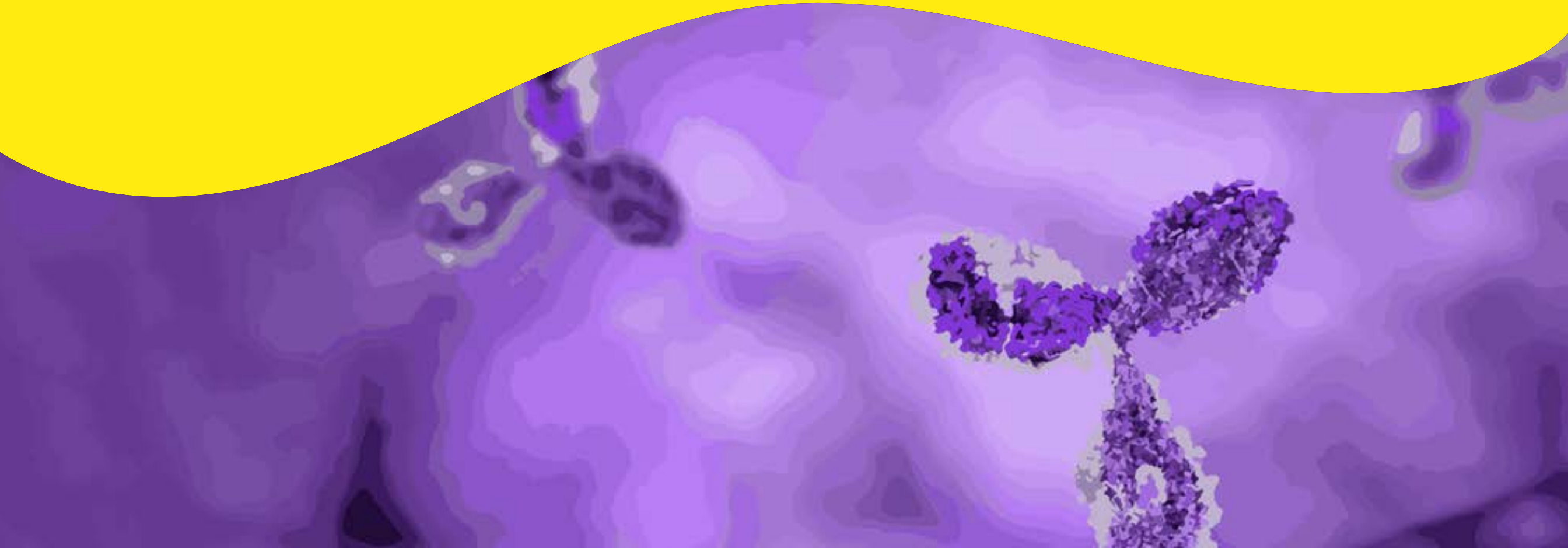


revvity

Autoimmunity

Guide to autoimmune diseases and therapeutic strategies



PURPOSE AND SCOPE

What to expect?

Welcome to this Autoimmune Diseases Booklet, a document to help scientists and researchers appreciate and navigate the diversity of the cell types and molecular pathways associated with autoimmunity, and the dysregulated immune responses associated with autoimmune diseases. We hope the visuals provided in this document will shed light on and clarify these otherwise complex mechanisms.

Among the over 80 autoimmune diseases reported so far, this document focuses on Rheumatoid Arthritis, Systemic Lupus erythematosus, Inflammatory Bowel Disease, Multiple Sclerosis, and Type I Diabetes.

You will find three separate sections:

The first groups the main facts understood about the pathogenesis underlying each of the autoimmune diseases listed above, with a short description of its associated symptoms and complications, as well as risk factors and key cellular and molecular involvements.

The second section dives more deeply into the cellular and molecular mechanisms generally associated with the diseases.

The last section gives an overview of current treatments used for autoimmune diseases, as well as some future directions.

The collection of cellular and molecular basics presented in this document was prepared based on authentic and highly regarded articles and journals. The numbers in brackets indicate the references used, and all pathways have been curated for scientific knowledge and accuracy by Revvity's scientific team.

Why this guide?

This new guide continues Revvity's tradition of providing collections of specialized documents dedicated to different therapeutic areas such as immunology, neurosciences, diabetes, and NAFLD. We also give you more directly practical guides covering the expertise of Revvity's scientists in assay development and performance.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
 - Systemic lupus erythematosus
 - Inflammatory bowel disease
 - Multiple sclerosis
 - Type 1 diabetes
 - Common antigens and autoantibodies associated with AI
-

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
 - Innate immunity components
-

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
 - Immunosuppressive drugs
 - Promising new therapeutic areas
-

Bibliography



AUTOIMMUNE DISEASES AT A GLANCE

Autoimmune diseases

Like an army of experienced and disciplined soldiers, a healthy immune system fights pathogens or abnormal cells to protect the body against diseases. But if the immune system dysfunctions, it may mistakenly mount an attack against healthy cells, tissues, and ultimately organs. These autoimmune diseases can affect any part of the body, weakening organism function and even becoming life-threatening.

AIDs mainly affect women, as 80% of the patients are women, and it has been calculated that there are around 300 million patients in the world. AIDs can be triggered by genetic and/or environmental factors, so the causes vary widely. In this booklet, we focus on the main autoimmune diseases: Rheumatoid Arthritis, Inflammatory Bowel Diseases, Atherosclerosis, Lupus and Type 1 diabetes.

To date more than 80 autoimmune diseases have been reported and classified in two categories: multisystem and organ specific autoimmune diseases. Multiple Autoimmune Syndrome (MAS) refers to the combination of several autoimmune disorders. It is known that 25% of patients suffering from one autoimmune disease tend to develop additional autoimmune diseases. For example, Rheumatoid Arthritis, Systemic Lupus erythematosus, Systemic Sclerosis, Sjogren’s syndrome, or spondyloarthritides are multisystem autoimmune diseases. Organ-specific autoimmune diseases refer to one organ or tissue preferentially attacked by the patient’s immune system. For instance, Type 1 diabetes patients suffer from pancreatic beta-cell destruction, and Vitiligo affects skin pigmentation, whereas Multiple Sclerosis affects the brain and spinal cord.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

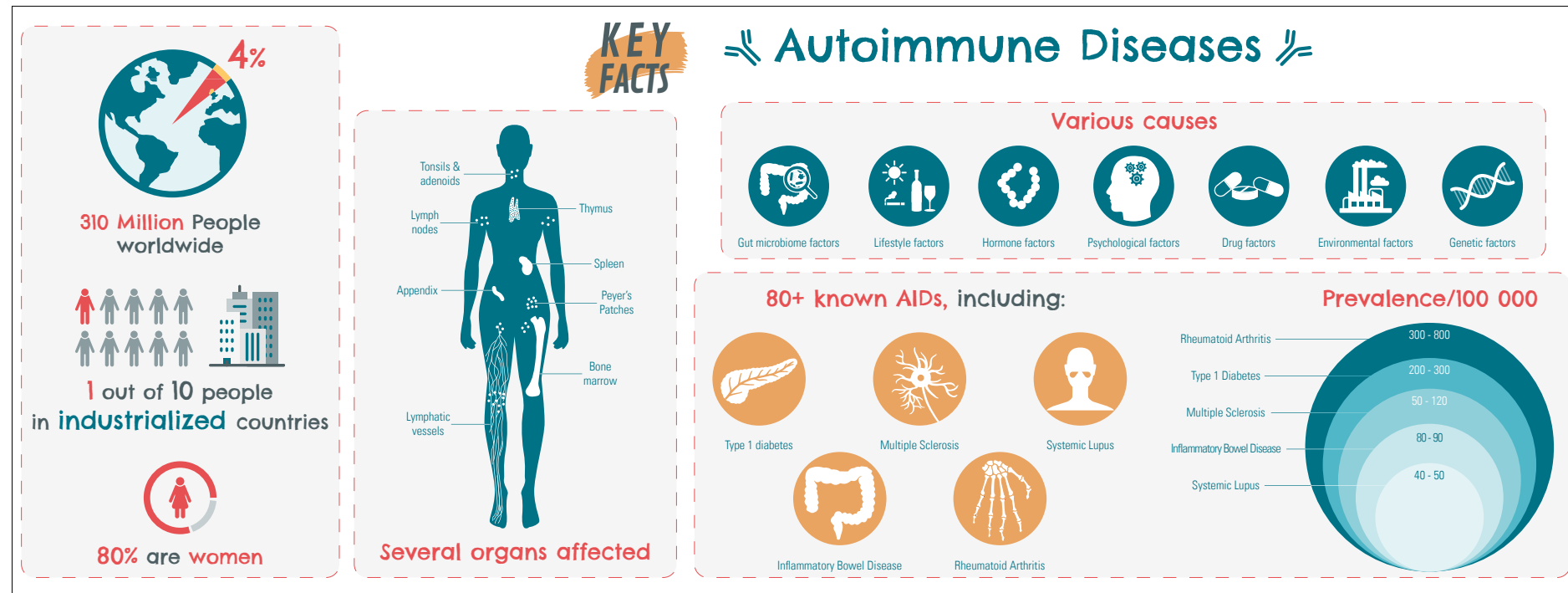
Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Rheumatoid arthritis

RA is the most common chronic inflammatory disease. This autoimmune disease affects around 1% of the population in the world, and is more prevalent in females than males (1).

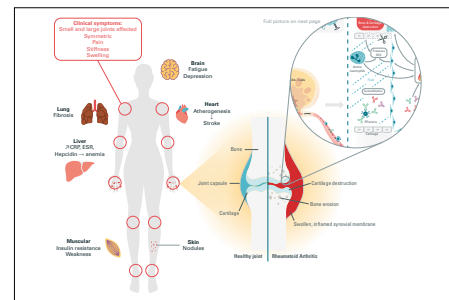
RA is a systemic and symmetric inflammatory disease associated with joint redness, hotness, stiffness, swelling and pain. Affected joints are primarily located at wrists, fingers, toes, and knees. In normal conditions, the synovium membrane surrounding the joints produces the synovial fluid which protects the joint. However, in RA inflammation of the synovium leads to an excess of synovial fluid. RA synovitis leads to cartilage damage, bone erosions, and eventually joint destruction.

Environmental and genetic factors play important roles in the disease onset. Among environmental factors, low socioeconomic conditions, diet and gastrointestinal microbiome, or smoking are the most recognized (2). Several genes have been associated with RA predisposition, such as the HLA-DRB1b gene involved in antigen presentation (3), the PTPN22 gene encoding for a protein tyrosine phosphatase involved in the negative control of T-cell receptor signal transduction, the TRAF1 gene involved in the negative regulation of signaling via TNF receptors 1 and 2 (4) and STAT4 implicated in the signaling of various cytokines like IL-12, IL-23, and type I interferons (5). In addition, epigenetic regulation of gene expression, including DNA methylation, histone phosphorylation, acetylation, has been postulated to play roles in RA pathology (6).

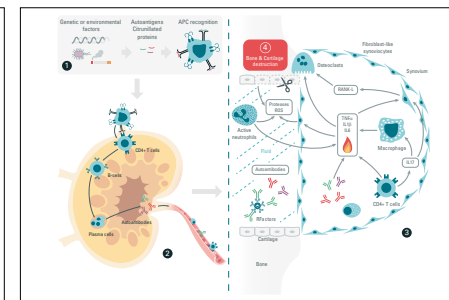
At the molecular level, genetic or environmental factors lead to the production of RA associated autoantigens, such as citrullinated proteins (7). Citrullination is a post-translational modification that converts arginine into citrulline. Citrullinated proteins, such as fibrin, fibrinogen, or vimentin, are then considered to be foreign antigens and are recognized by Antigen-Presenting Cells. Autoantigen loaded-APCs migrate into the lymph nodes where they activate CD4+ helper T-Cells. In turn, these helper T-cells stimulate B-cell differentiation into plasma cells which produce autoantibodies. The two most important autoantibodies associated with RA are Rheumatoid Factor (RF) which are essentially IgM and bind to the Fc portion of IgG, and anti-citrullinated protein antibodies (ACPA) (8). Finally, CD4+ Helper T cells, plasma cells, RF, and ACPA auto antibodies migrate through the blood stream and infiltrate back into the joint.

At the joint, fibroblast-like synoviocytes constituting the synovium membrane, as well as macrophages, play a central role in the inflammation of the synovium called synovitis. Macrophages produce proinflammatory cytokines such as TNF α , IL1b, and IL6, which lead to inflammation. In addition, cytokine-activated synoviocytes secrete RANK-L and metalloproteinases. Together, cytokines and RANK-L promote the activation of the osteoclasts leading to bone erosion, whereas metalloproteinases are essentially responsible for cartilage degradation. Furthermore, infiltrated activated CD4+ T-cells secrete IL17 and RANK-L contributing to macrophage, synoviocyte, and osteoclast activation. Accentuation of bone and cartilage degradation also results from neutrophils producing proteases and reactive oxygen species (9). Moreover, synovitis associated Plasma cells participate in inflammation through cytokine and autoantibody production. Here, immune complex deposition of RF and ACPA autoantibodies with complement proteins leads to a severe inflammation. Finally, increased angiogenesis and vascular permeability reinforce immune cell infiltration into the joint.

In addition to attacks on the joints, the functionality of several organs is impaired in RA due to elevated concentrations of TNF α , IL1b, and IL6 in the circulation. The extra-articular manifestations reside i) in the liver, producing c-reactive protein and hepcidin-induced anemia, ii) in the skin, by nodule formations, iii) in the heart, by atheroma plaque formation, iv) in the central nervous system, by fatigue and depression, v) in the muscles, by insulin resistance, and vi) in the lungs, by fibrosis.



[Click to Enlarge](#)



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

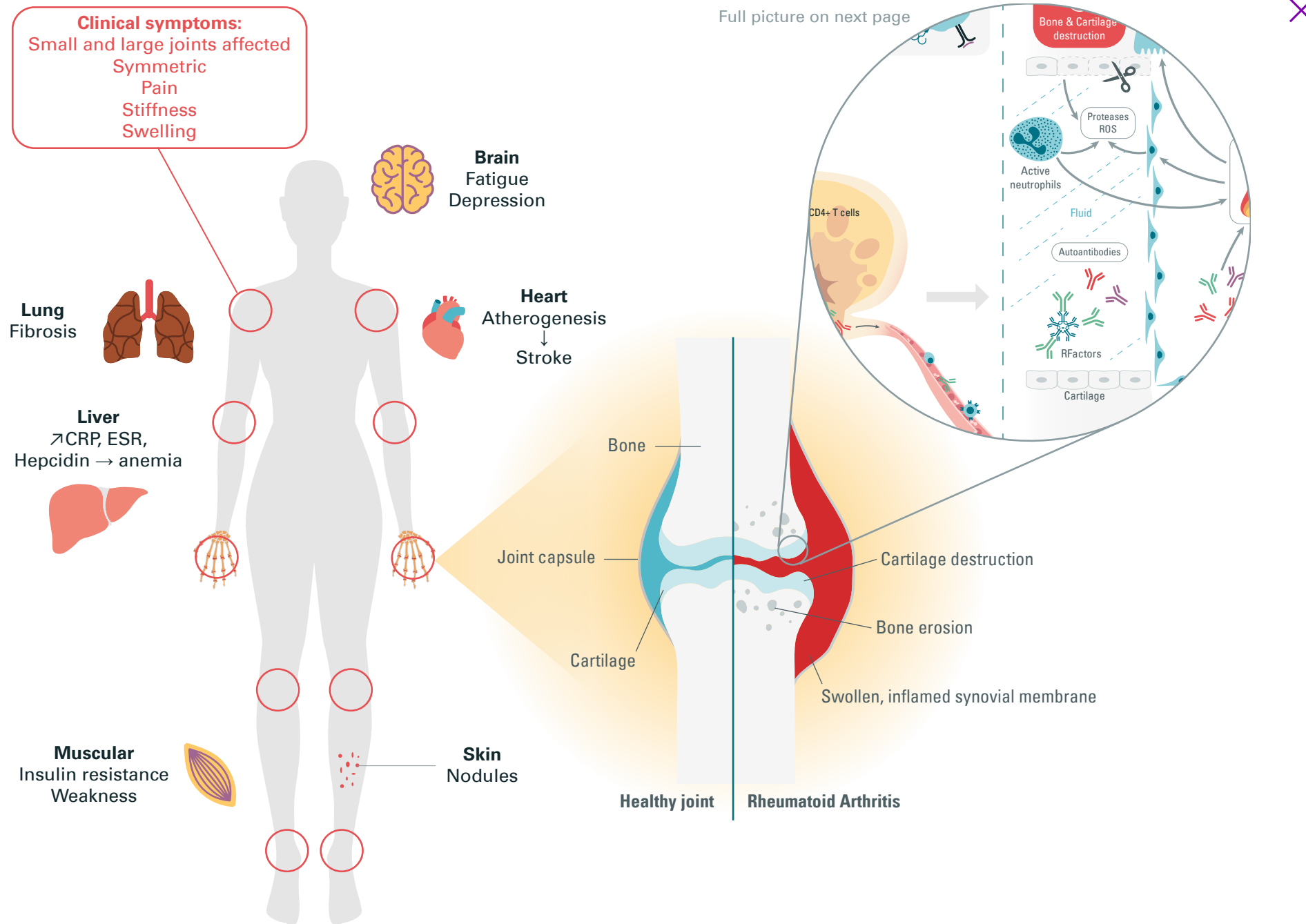
Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



Full picture on next page



TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

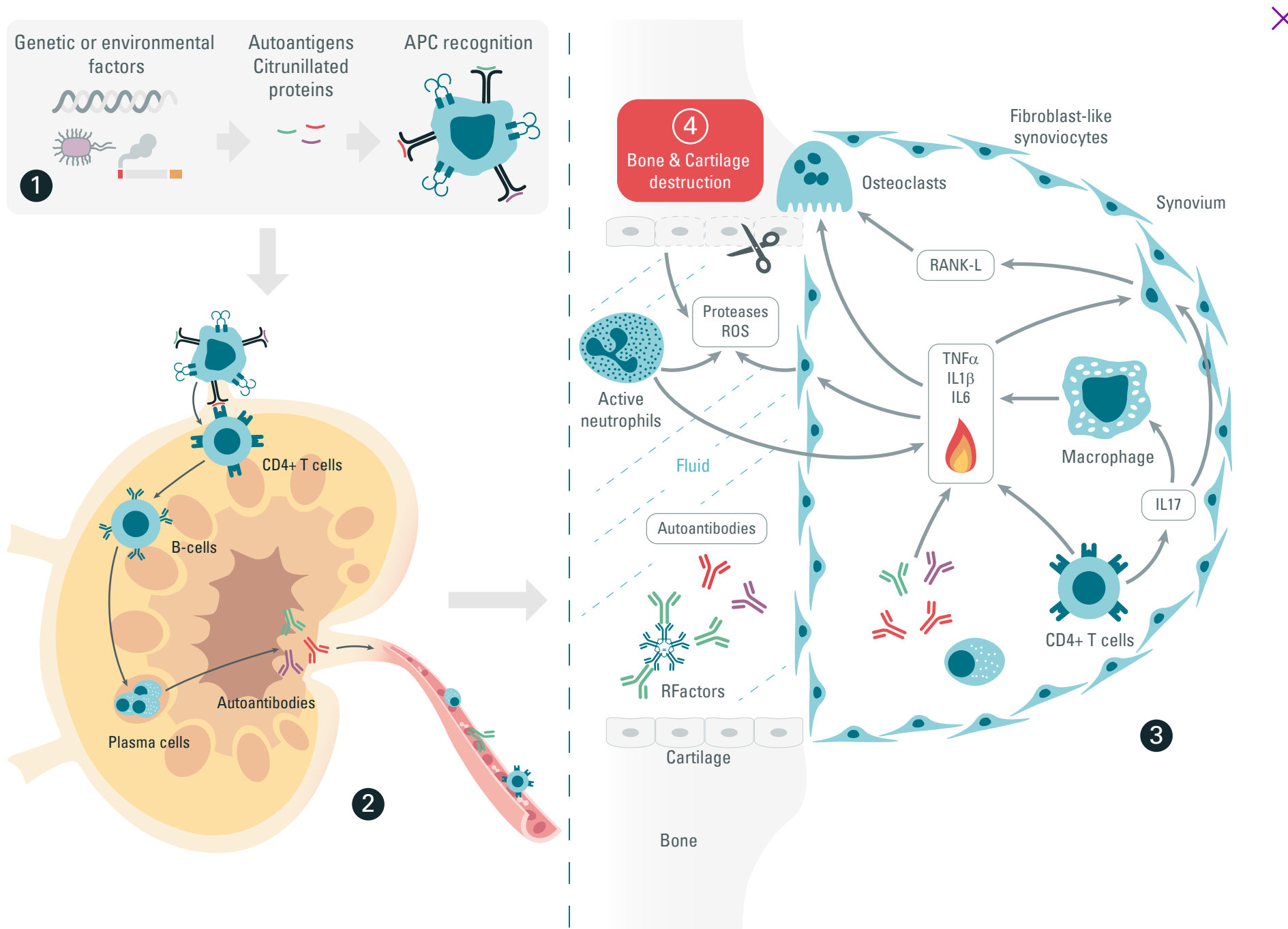


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Systemic lupus erythematosus

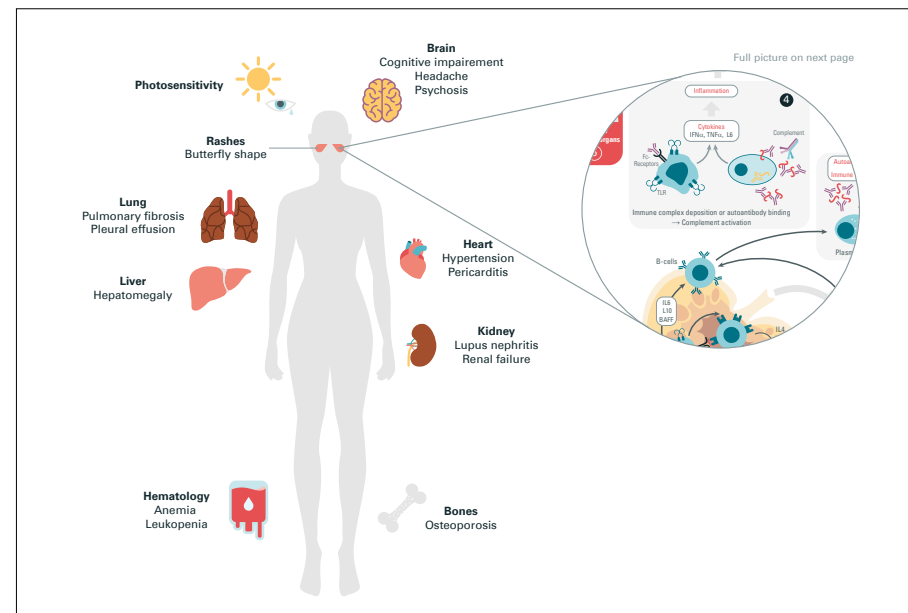
Systemic Lupus Erythematosus, or SLE, is an autoimmune disease mediated by autoantibodies and immune complexes. SLE is a chronic multisystem disorder mainly affecting women between 15 and 40 years of age (10). Symptoms associated with SLE are fatigue, weight loss, and fever. UV sensitive butterfly rash over the nasal bridge occurs in 50% of SLE patients and represents the most common sign of SLE. Neurological affection is common and includes cognitive impairment and psychosis, as well peripheral neuropathies. Pulmonary fibrosis, vasculitis, and pleuritis are also experienced in some SLE cases, as well as heart manifestations causing chest pain and breathing issues. Lupus nephritis and ultimately renal failure result in hypertension which is a common complication of SLE. Other symptoms such as hepatomegaly, abdominal pain, nausea, chronic anemia, polyarticular arthritis, or osteoporosis are also associated with it.

SLE results from interplay between genetic, immunological, and hormonal factors, as well as environmental factors. For example, deficiencies in early complement proteins such as C1q and C4 are associated with SLE, as well as SLE-associated gene variants (e.g. IRFs, STAT4, or BLK) (11) (12). Mutations in the intracellular DNA exonuclease, TREX1, in SAMHD1 or ADAR1 genes have also been associated with lupus (13), leading to the accumulation of nucleic acids and the subsequent activation of the innate immune system. SLE has also been related to elevated Estrogen during pregnancy (14).

Environmental factors may act as initial triggers of SLE. These include UV light, viral infections such as Epstein-Barr Virus, and smoking. Environmental factors induce DNA damage and ultimately cell death through apoptosis or necrosis (15). Along with "neutrophil extracellular traps", also called NETosis, nuclear as well as cytoplasmic and membrane components are released and exposed to the innate cells (16). Immature antigen presenting cells sense cellular material through Toll Like Receptors (TLR) or nucleic acids sensors, such as cGAs or RIG-I proteins (17) (18), which in

turn induce high levels of Type-I IFNs. Nuclear antigens include Histone core, dsDNA, and ribonucleoprotein complexes including SM, VI RNP, and Ro/SS-A. Cytoplasmic components also include Ro, whereas cell membrane antigens include cardiolipin found in mitochondria, platelets, or red blood cell membrane components (19).

Regardless of the antigens that are picked up by the APC, antigens are presented to naïve CD4+ T helper cells in the lymph nodes. Activated helper T-cells, mainly Th2, Th17, and follicular helper T cells, promote B cell activation, proliferation, and differentiation into plasma cells (20) (21). Then the plasma cells produce autoantibodies, such as anti-dsDNA antibodies and anti-histone antibodies, also known as ANA or Anti-Nuclear antibodies. So far more than 100 autoantibodies have been reported in SLE patients (22). These autoantibodies form immune complexes that can target virtually any organ.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

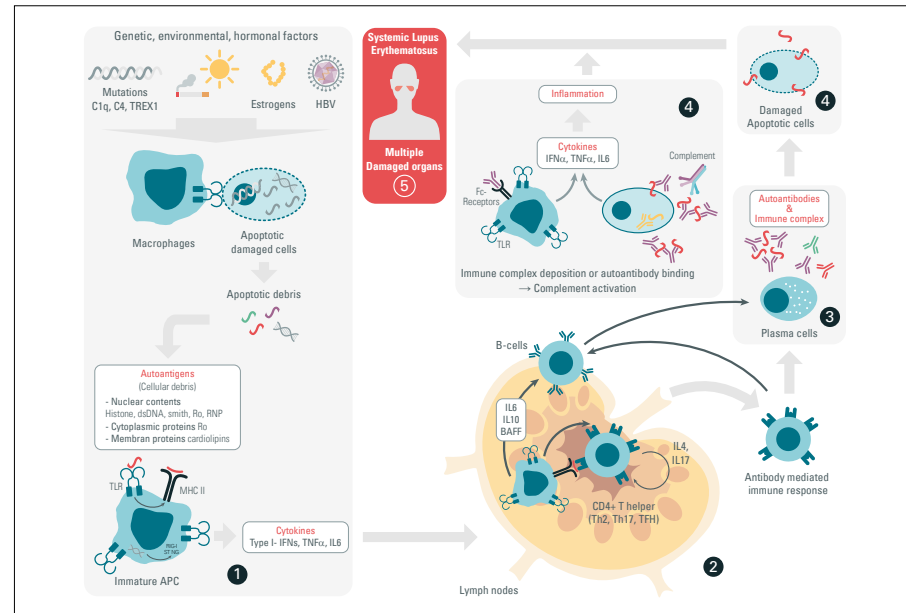
Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Systemic lupus erythematosus continued

Several mechanisms participate in the inflammation process which ultimately leads to tissue injury. Immune complex formation and deposition in organs results in complement activation and inflammatory response. The immune complexes can also bind to the Fc receptors of immune cells which will trigger the release of type I IFNs and proinflammatory cytokines such as TNF α or IL6, thereby increasing inflammation. The autoantibodies may bind on antigens present on the non-hematopoietic cell surface, resulting in complement activation and further cytokine releases. Finally, the sensitized immune cells detect antigens through innate immune sensors (e.g. TLR) and release proinflammatory cytokines which exacerbate the inflammation. Thus, amplified inflammation causes further organ injury, producing more damaged cells, and leading to an endless cycle.

SLE is an on-going chronic disease associated with flare-ups which reflect an immunological B cell memory and occur in response to a new exposure to antigens. Any precipitating factors causing cell apoptosis, such as sun exposure, infections, stress, or pregnancy, can trigger flare-ups and exposure to the nuclear antigens. SLE patients may have increased levels of B-cell activating factor BAFF involved in B cell survival and maturation, and antibody production by plasma cells, thus contributing to the persistence of autoreactive B cells and immunological B cell memory (23).



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

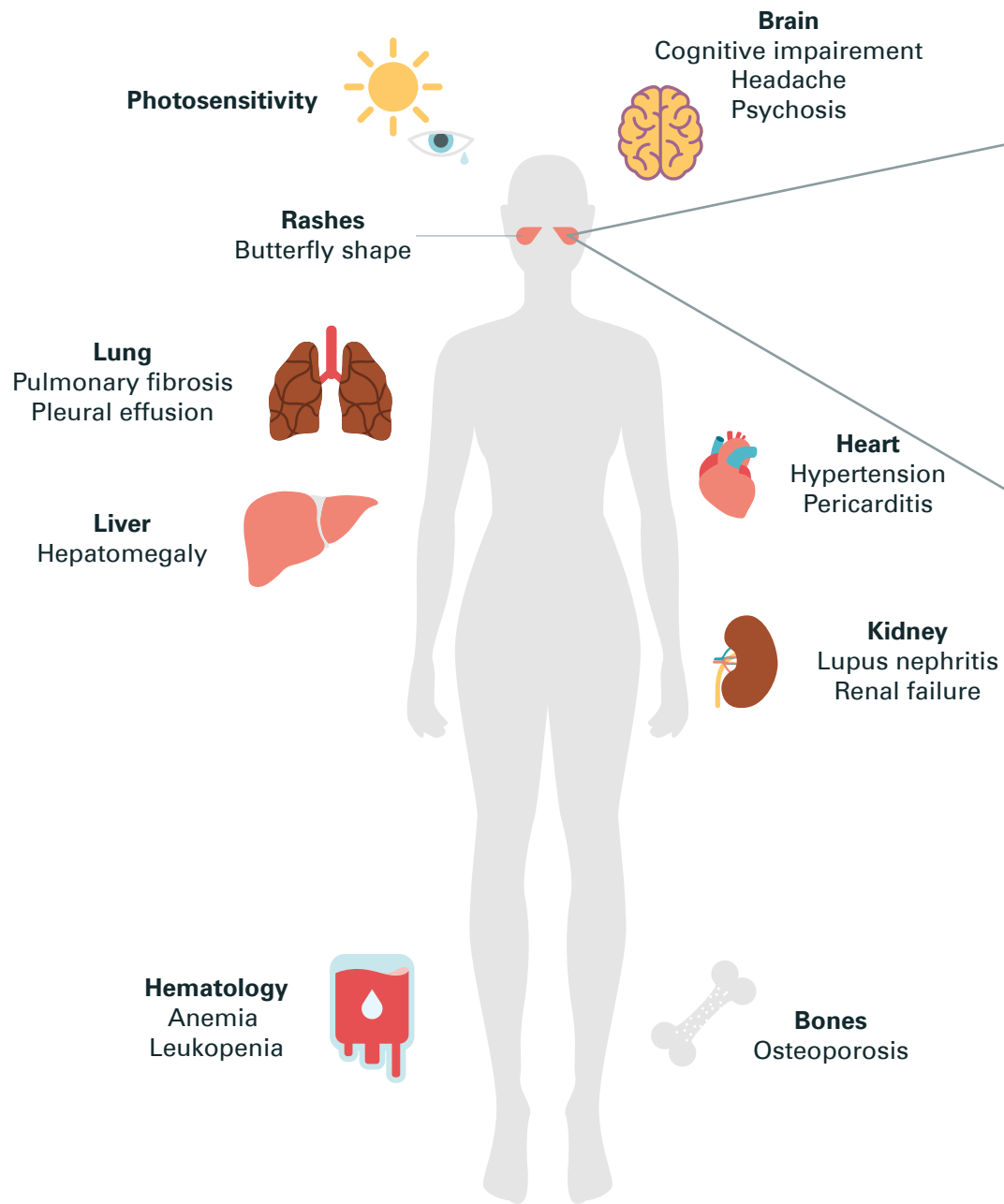
Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



Full picture on next page

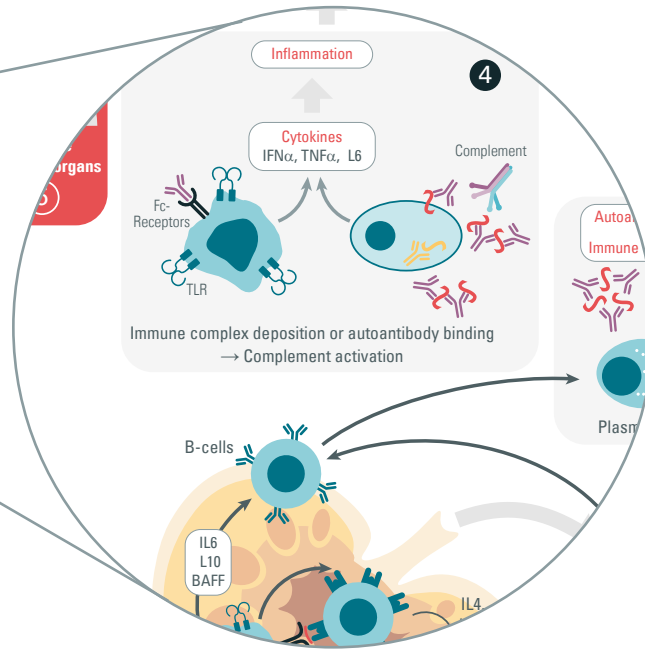


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

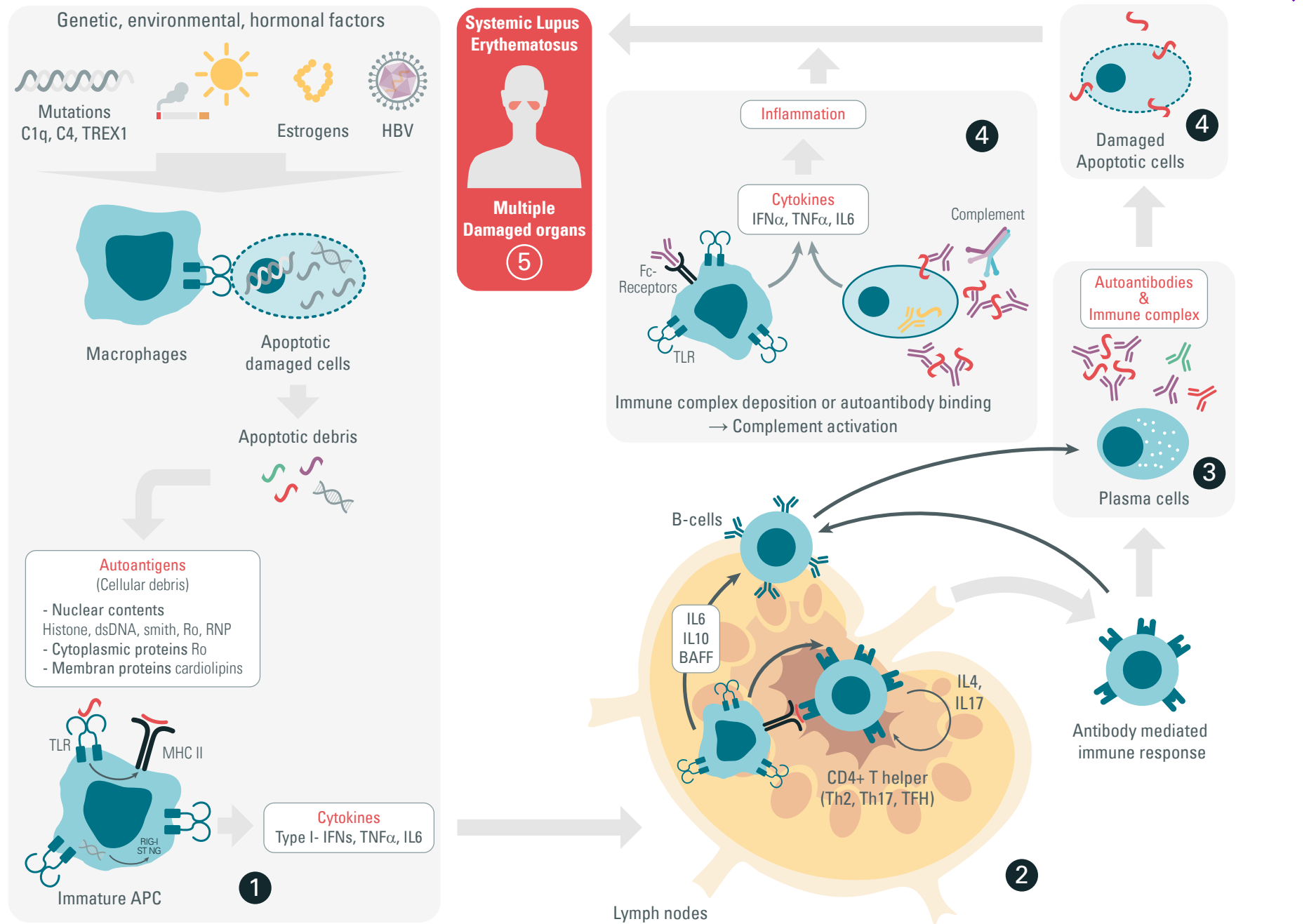


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

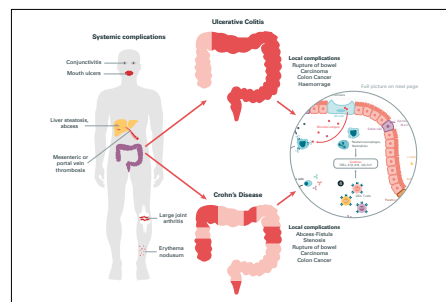
Inflammatory bowel disease

Inflammatory bowel disease (IBD) is divided into two related diseases: Crohn's Disease (CD) and ulcerative colitis (UC), and often appears in young adults between 15 and 40 years of age. UC is a continuous superficial inflammation that spreads from the distal to the proximal colon and is mainly localized in the rectal and sigmoid colon. CD is a discontinuous inflammation affecting the different cellular layers of the gastrointestinal tract, mainly the ileocecal area, the small intestine and colon. Symptoms associated with IBD can include diarrhea, rectal bleeding, abdominal pain, weight loss, and cramps. Complications associated with UC include severe bleeding, polyps, and rupture of the bowel, as well as colon cancer. CD complications include stenosis, abscess, fistula, fibrosis, and colon cancer, as well as perforation. IBD is associated with several extraintestinal manifestations, involving liver steatosis, thrombosis, joint arthritis, conjunctivitis, or mouth ulcers.

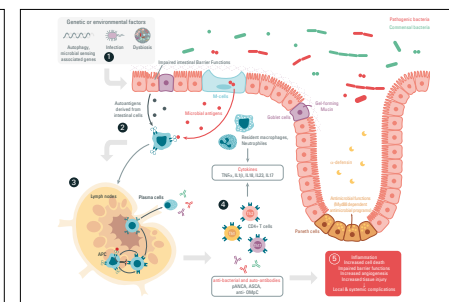
Combinations of environmental and genetic factors contribute to IBD development. For example, diet, stress, virus infections, non-steroidal anti-inflammatory drugs, or antibiotics, and smoking can promote IBD, as well as some gene variants encoding cytokines or their downstream signaling mediators (24), autophagy (e.g. ATG16L) (25), and microbial sensing (e.g. NOD2 gene) (26) (27). Dysbiosis, which refers to microbiota perturbations, has also been associated with IBD (28). In healthy patients, commensal microorganisms (e.g. *B.fragilis*) ensure immune protection by promoting the generation of peripheral T-reg cells and the subsequent production of the anti-inflammatory cytokine IL10. IBD patients show a high abundance of *Escherichia Coli* and other bacteria that promote inflammatory responses. In addition, in CD patients the mutation of the intracellular bacterial sensor NOD2 gene is associated with inflammatory gut microbiota.

Impairment of the epithelial barrier function and dysbiosis are key events in IBD initiation. Microbial peptides, which are normally translocated through M cells, pass through the lumen due to intestinal barrier impairment. Microbial and intestinal epithelial cell (IECs)-derived antigens activate APC, such as macrophages and dendritic cells, which promote CD4+ T helper cell activation. IECs, activated macrophages, and CD4+ T lymphocytes secrete proinflammatory cytokines such as

TNF α , IL1b, IL18, IL17, and IL23, as well as IFN γ , IL6, and IL33, resulting in chronic inflammation (29) (30) (31). In addition, antigens derived from IECs lead to B cell activation, proliferation, and differentiation into plasma cells. Autoantibodies secreted by plasma cells (Anti-Neutrophil Cytoplasmic Antibody :ANCA), along with the anti-bacterial antibodies Anti-Saccharomyces Cerevisiae Antibody (ASCA) or anti-Outer-membrane protein of *Escherichia Coli* (anti-OMPc) participate in the inflammatory process through the formation of immune complexes and complement activation, or the activation of macrophages through Toll-like receptors (32). Taken together, the activation of the immune system results in local and systemic complications where TNF α plays pleiotropic roles in IBD pathogenesis (33). TNF α stimulates angiogenesis, increasing the presence of immune cells locally. TNF α induces the necrosis of Paneth cells, as well as Goblet cells, halting antimicrobial functions and mucus production respectively. TNF α induces Intestinal Epithelial Cell death and impairs intestinal barrier functions. TNF α increases immune system activity, thus causing further intestinal damage. Finally, TNF α stimulates myofibroblasts to release proteases which further destroy the intestinal epithelium.



[Click to Enlarge](#)



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

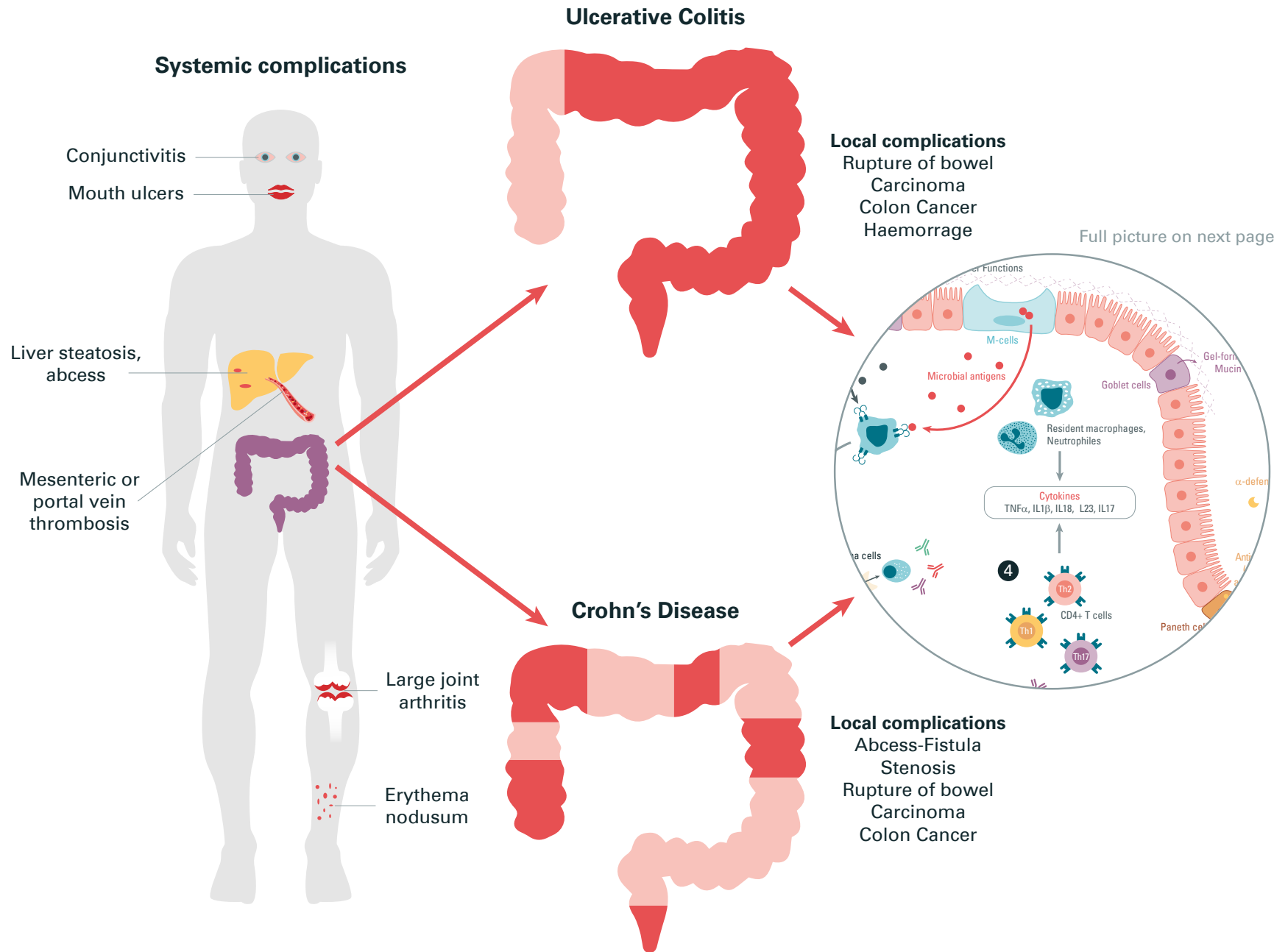


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

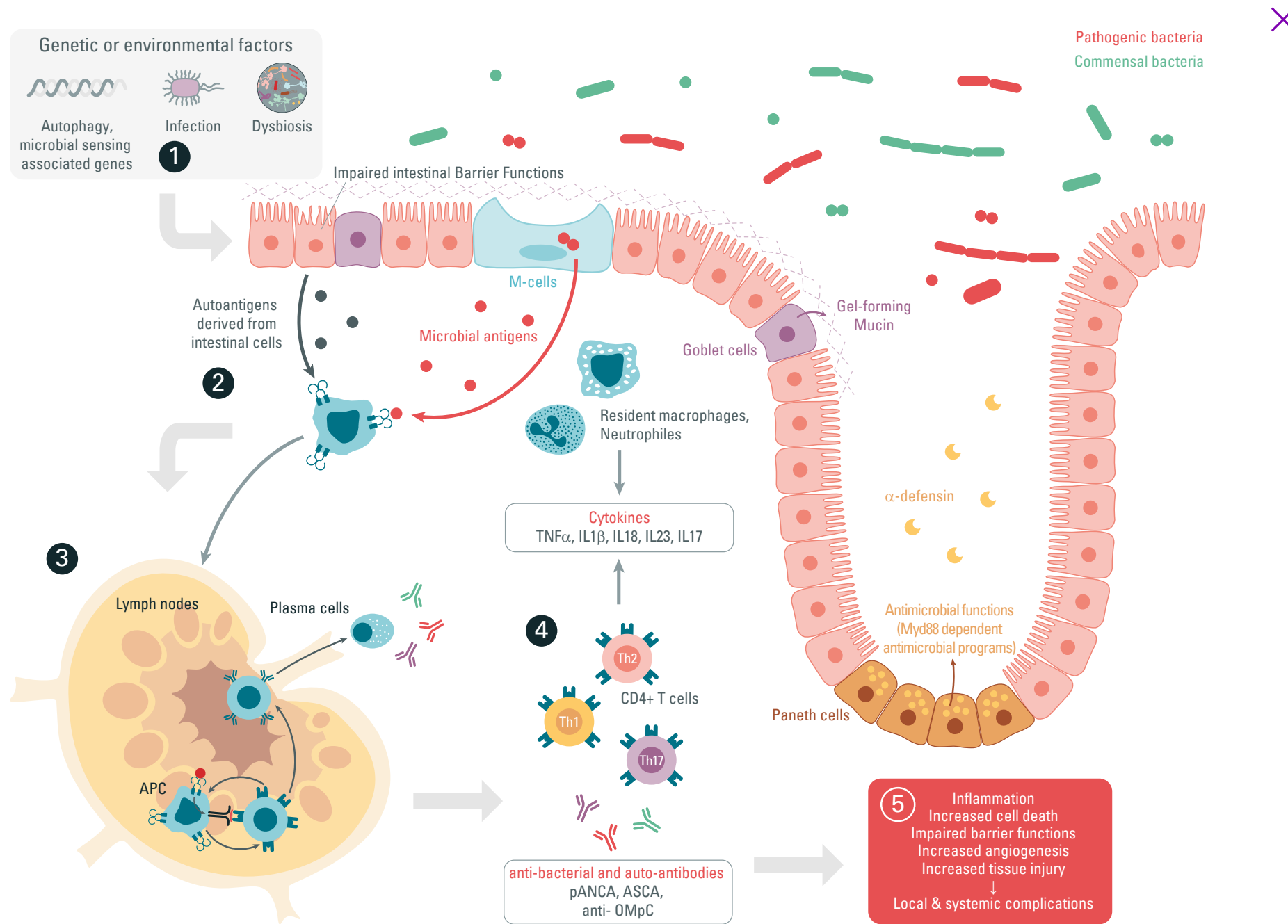


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Multiple sclerosis

MS is a chronic organ-specific autoimmune disease targeting the Central Nervous System. MS is associated with neuroinflammation resulting in a progressive demyelination process and neuronal damage. Typical symptoms include temporary visual loss due to optic neuritis, motor and sensorial troubles, pain, fatigue, depression, neurocognitive troubles, but also impairment of bladder, bowel, and sexual functions (34). Relapsing-remitting and secondary progressive MS (RRMS, SPMS) are the two most frequent forms of MS, predominantly affecting women between 20 and 40 years of age. RRMS is marked by sequential relapses or exacerbations of symptoms followed by periods of remission, when symptoms improve or disappear. In patients suffering from SPMS, the course of the disease progressively worsens with or without periods of remission. The Primary progressive MS (PPMS) form is more resistant to treatment and symptoms become steadily worse without remissions. Progressive-relapsing MS (PRMS) is less common, progressive, and associated with flare-ups of symptoms with no periods of remission.

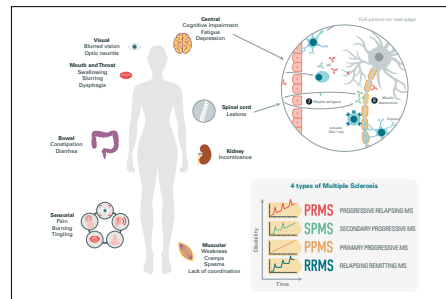
Despite decades of investigations, the initial trigger of MS still remains elusive. Nevertheless, the interplay between genetic, epigenetic, and environmental factors is likely to play a role in the onset and the progression of MS (35). Whereas the involvement of epigenetically regulated allele variant HLA DRB*1501 has recently been found in MS pathogenesis (36) (37), the HLA-A2 allele seems to be protective. Viral infections, particularly with herpes virus, low Vitamin D due to reduced sun exposure, diet and obesity, gut microbiota, air pollution, but also smoking have been identified as contributing to MS (38) (39).

The perturbation of peripheral immune responses, as well as brain cells such as microglia or astrocytes, convey an autoimmune process to the CNS.

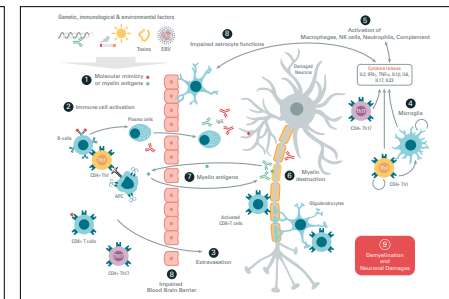
Whether the antigens inducing immune activation are molecular mimicry (viral or bacterial peptides that share homology with brain proteins such as myelin) or myelin antigens (40), activated CD4+ Th1, CD4+ Th17, and CD8+ peripheral T cells infiltrate the CNS where they secrete pro-inflammatory cytokines such as IL2, IFN γ , TNF α , or IL6. CD8+ cytotoxic T cells directly attack the oligodendrocytes and the myelin, promoting their destruction (41). Activated microglia, macrophages, NK cells, or neutrophils participate in pro-inflammatory cytokine secretion and thereby

contribute to neuronal damage and a demyelination process. In addition, activated B cells secrete autoantibodies which amplify the CNS attack, through complement activation or macrophages inducing neuron damage (41). Of note, even though 90% of the patients have elevated levels of CSF IgG, only a few specific autoantibodies have been attributed to MS, with Anti-myelin oligodendrocyte glycoprotein (MOG) being one them (42) (43) (44). Furthermore, a continuous Myelin antigen release induces myelin autoantibody production and further amplifies the demyelinating circle. Finally, impaired astrocytes and blood brain barrier functions contribute to disease progression.

For more information regarding neuroinflammation and neurodegenerative diseases, please read our Neuroscience guide.



[Click to Enlarge](#)



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

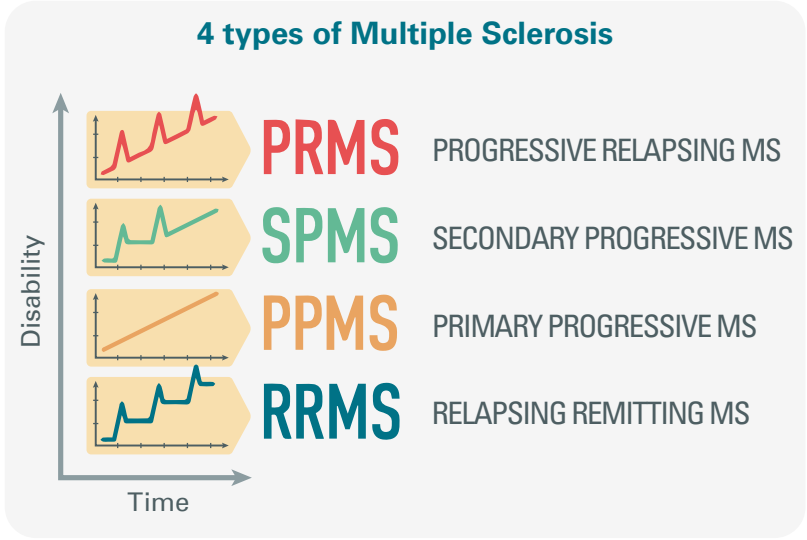
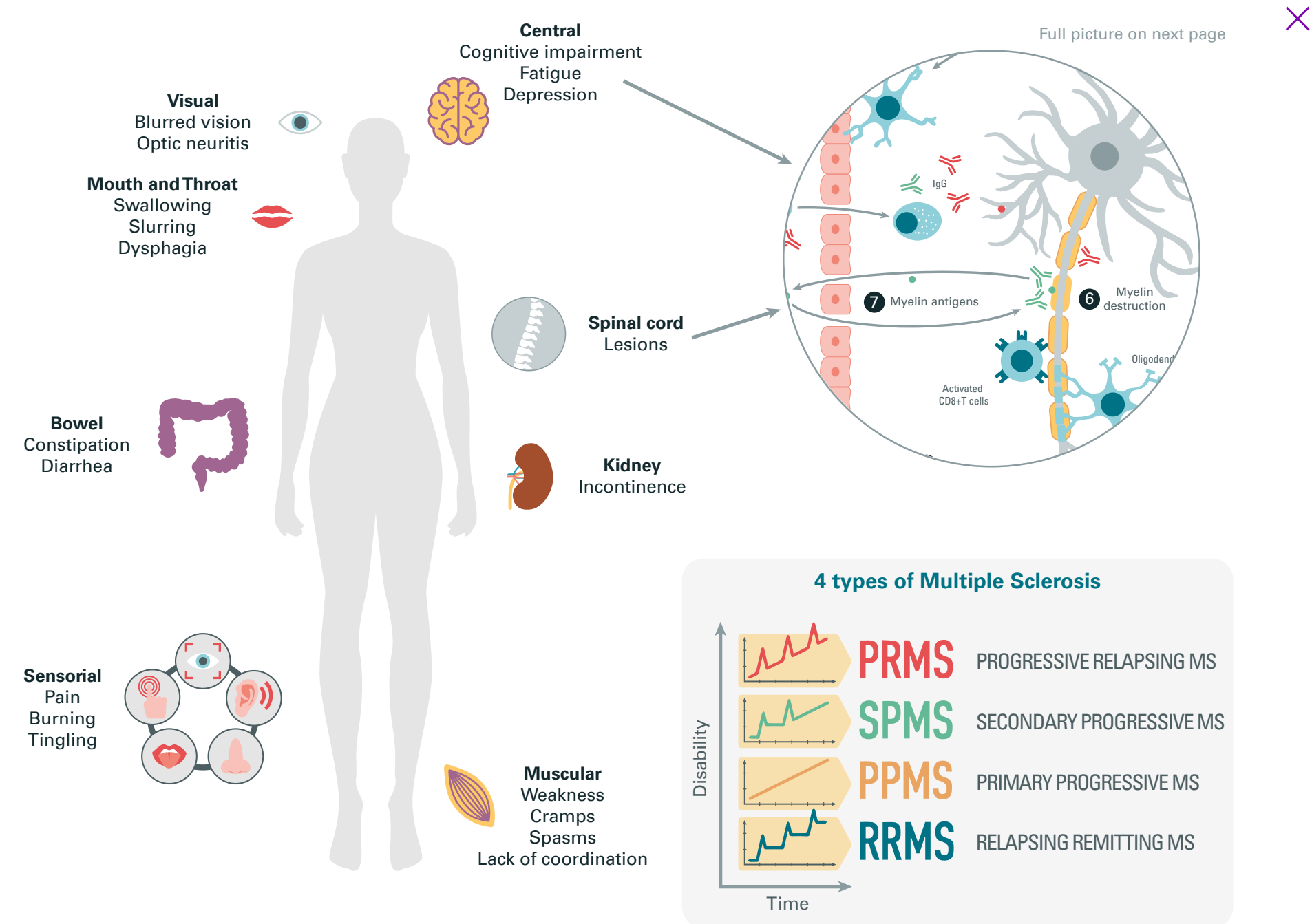


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



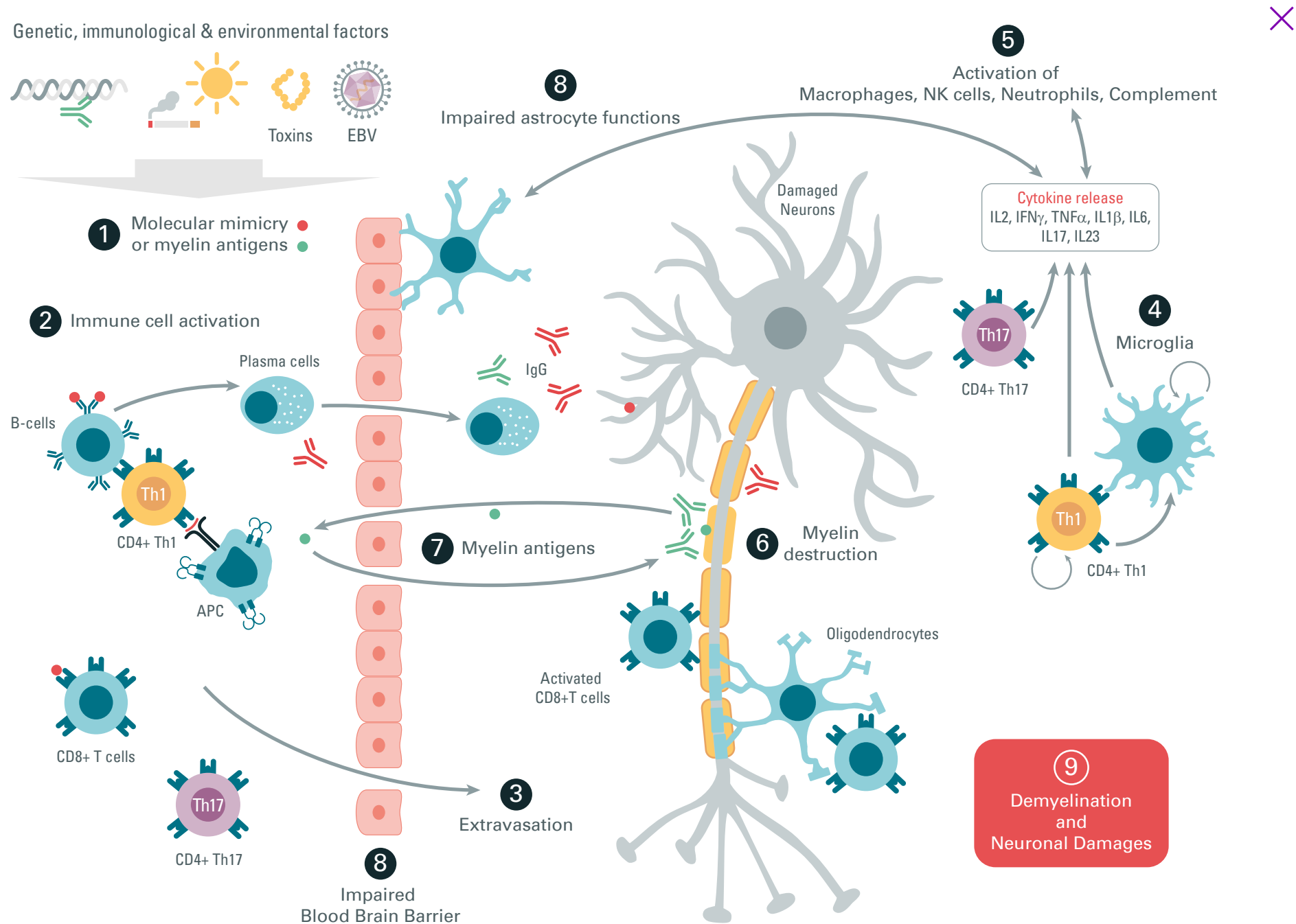


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Type 1 diabetes

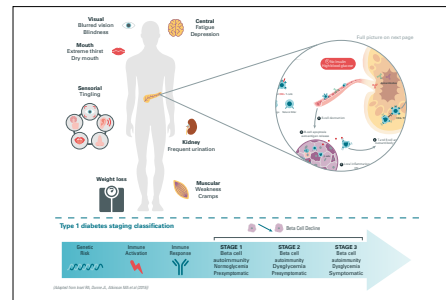
Type 1 diabetes is a chronic autoimmune disease, less common than its sibling Type 2 diabetes. Autoimmune diabetes represents 10% of diabetes patients and is the most common form affecting children under 15 years of age. Mellitus diabetes is characterized by high blood glucose levels due to pancreatic islet beta-cell destruction and insulin deficiency.

The pathogenesis of autoimmune type 1 diabetes is divided into three stages (45). Stage 1 is presymptomatic, the blood glucose level being normal (normoglycemia), and relates to β -cell autoimmunity with the presence of islet autoantibodies. Stage 2 is still presymptomatic, but glycemia is perturbed, and finally stage 3 sees the onset of symptomatic disease. Typical symptoms of diabetes include frequent urination, increased thirst, hunger but weight loss, fatigue and irritability, but also blurred vision and skin infections. Left untreated, type 1 diabetes can lead to life-threatening complications through diabetic ketoacidosis (DKA) such as neuropathy, nephropathy, heart and blood vessel disease, and coma.

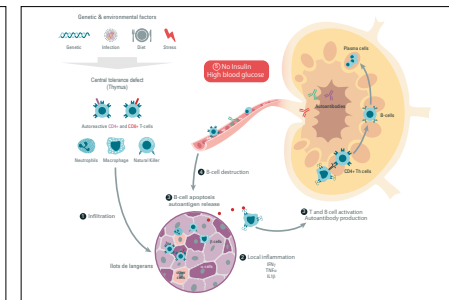
Even though the initial triggers of T1D pathogenesis remain to be elucidated, the interplay between genetic factors, mainly HLA alleles (46), central and peripheral tolerance, as well as environmental factors contribute to the disease onset. Autoreactive T cells, neutrophils, and NK cells as well as macrophages infiltrate the beta-pancreatic islets, leading to inflammatory cytokine production such as TNF α , IL1b, and IFN γ , and local inflammation. Pancreatic insulin-producing β cells undergo

apoptosis, liberating autoantigens which are then recognized by Antigen Presenting Cells and migrate to the lymph nodes. Finally, activated CD4+ and CD8+ T cells, as well as autoantibody-producing differentiated B cells leave the lymph node and migrate back to the pancreas, where they exert their cytotoxic activities, leading to β -cell destruction and T1D (47). Major islet autoantibodies target insulin (IAA), GADA, islet antigen-2 (IA-2A), or the zinc transporter 8 (ZnT8A) (48).

For more information regarding diabetes and associated signaling pathways, please read our Diabetes guide.



[Click to Enlarge](#)



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

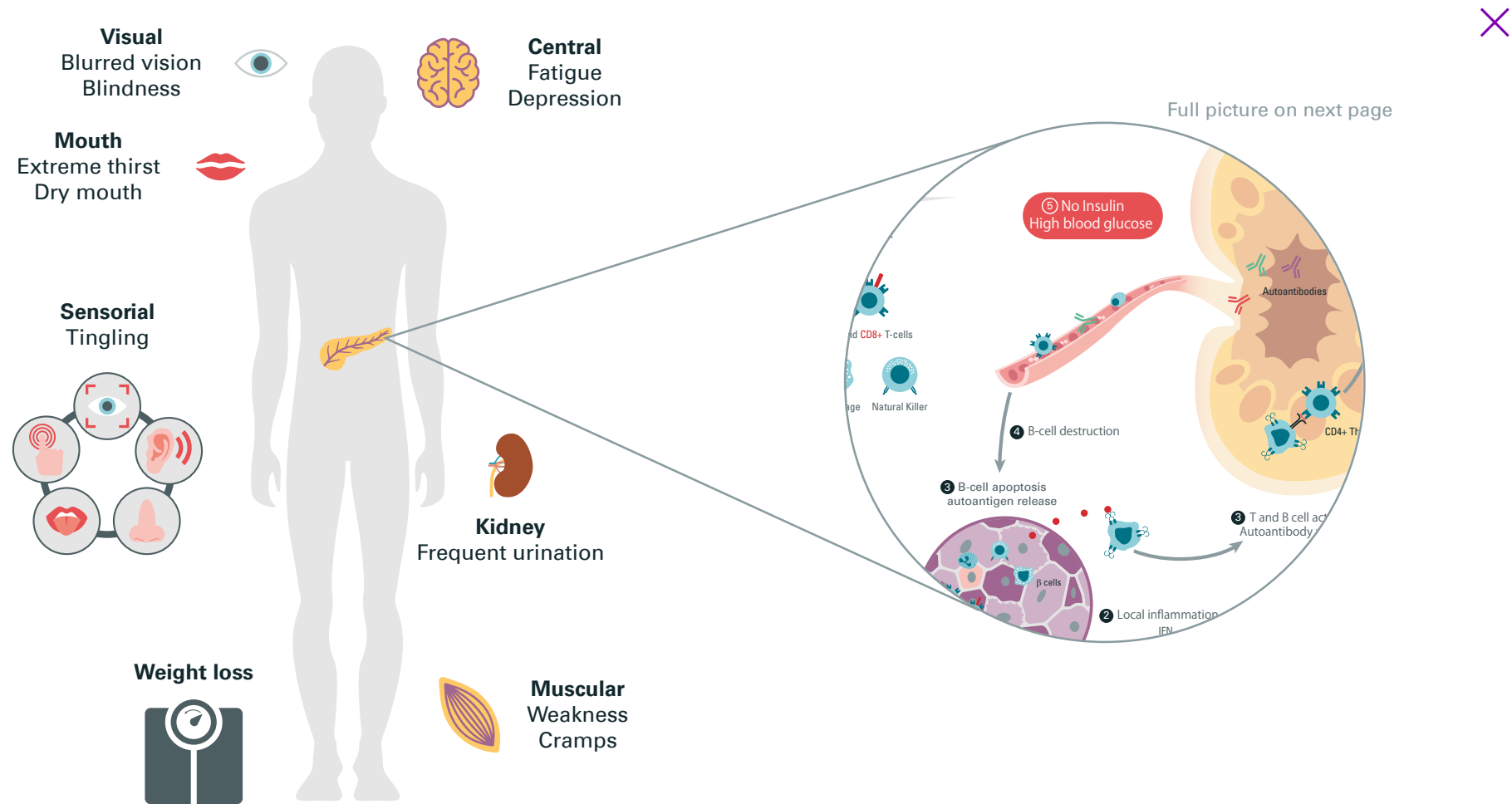
Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

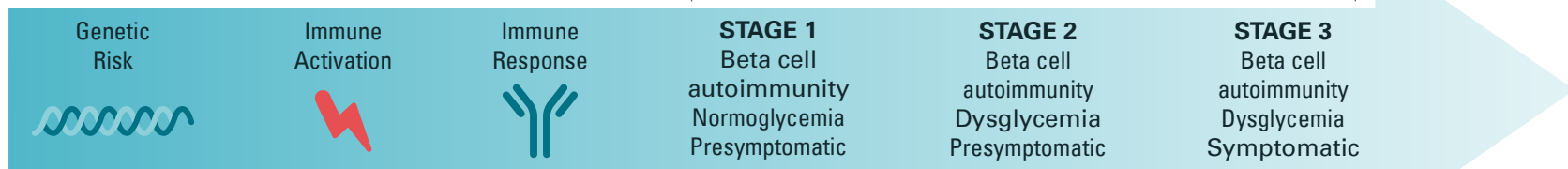
- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



Full picture on next page

Type 1 diabetes staging classification



(Adapted from Insel RA, Dunne JL, Atkinson MA et al (2015))

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

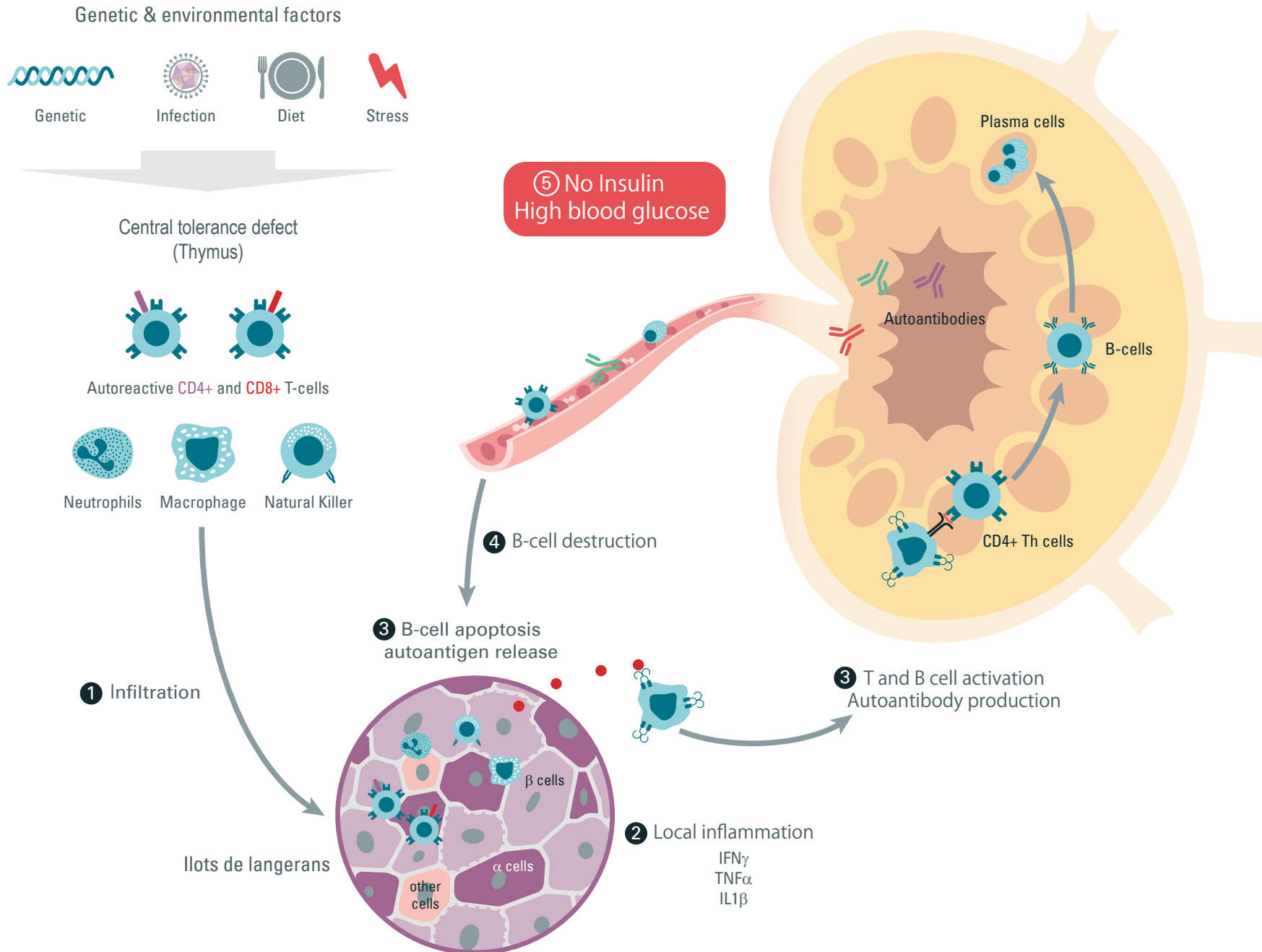


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

PATHOPHYSIOLOGY OF AUTOIMMUNE DISEASES

Common antigens and autoantibodies associated with AI

At some point, autoimmune diseases are elicited by antigens, whether self or microbial antigens, so-called molecular mimicry. These antigens can be captured by Antigen Presenting Cells, leading to the production of autoantibodies which in turn result in harmful effects in tissues.

The main autoantibodies associated with the previously described autoimmune diseases, as well as their targeted components, are listed in the following table:

AUTOIMMUNE DISEASES	AUTOANTIBODIES	TARGETED COMPONENTS
<i>Rheumatoid Arthritis</i>	Rheumatoid factor (RF) Anti Citrullinated Protein Antibodies (ACPA) anti-RA33	Fc portion of IgG. Fibrinogen, vimentin, filaggrin, CII, enolase. hnRNP
<i>Systemic Lupus Erythematosus</i>	Anti Nuclear antibody Anti-dsDNA Anti-Sm Anti-Ro/SSA Anti-La/SSB Anti-Cardiolipin	Nuclear components Double stranded DNA snRNPs complex hY-RNP complex Calmodulin-Binding Protein Mitochondrial protein
<i>Inflammatory Bowel Diseases</i>	perinuclear AntiNeutrophil Cytoplasmic Antibodies	Histone H1
<i>Multiple Sclerosis</i>	Anti-MOG Anti-AQP4	Myelin Oligodendrocyte Glycoprotein (MOG) Aquaporin-4 (water channel protein)
<i>Type 1 Diabetes</i>	Anti-IAA Anti-GADA Anti-IA-2A Anti-ZNT8A	Insulin Glutamic Acid Decarboxylase Islet antigen-2 Zinc transporter 8

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

CELLULAR AND MOLECULAR MECHANISMS OF AI DISEASES

Central and peripheral tolerance

In a healthy condition, our immune system does not recognize self-antigens and cannot mount an attack against our own tissues. This is known as “self-tolerance”, encompassing immune central and peripheral tolerance. The central tolerance mechanism occurs in lymphoid organs: the thymus for T-cells and the bone marrow for B-cells, where T or B lymphocytes that recognize self-antigens are eliminated or inactivated. Peripheral tolerance occurs in peripheral lymphoid organs, such as lymph nodes and spleen.

CENTRAL TOLERANCE

T-lymphocytes, produced in the bone marrow, are matured in the thymus, where different spatial and temporal steps orchestrate the differentiation and selection process (49) (50) (51) (52) (53).

In the capsular region, double negative T-cells (DN) express mainly TCRab but not the CD4 or CD8 co-receptors, and undergo TCR gene rearrangement, which provides the diversity of the TCR repertoire.

Then the DN cells move down to the cortex where they express both CD4 and CD8 co-receptors, now being called double positive T-cells (DP). Here DP interact and recognize a large panel of self-antigen peptides displayed on MHC-I or II expressed at the cell surface of thymic epithelial cells (TEC). This can be considered as the first degree of T-cell education and editing. Different selection procedures occur:

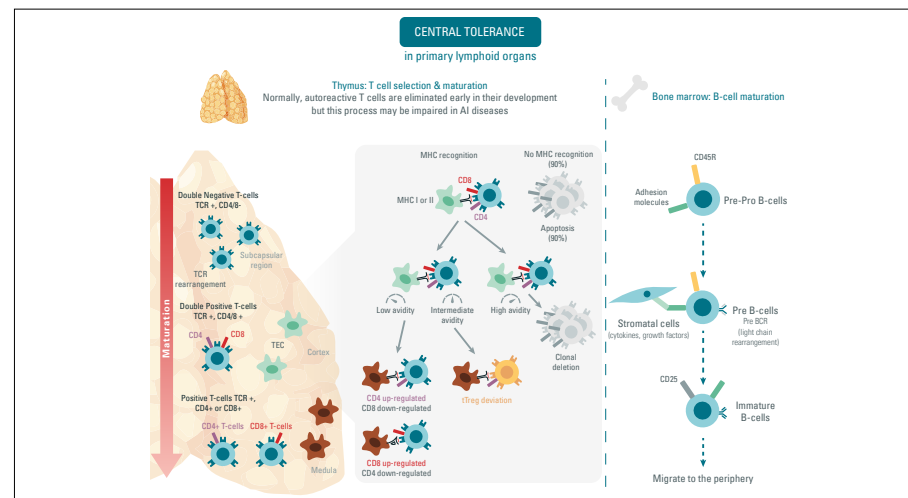
- A positive selection ensures T-cells strictly recognize peptide bound MHC; otherwise they undergo cell death, which occurs for 90% of T-cells.
- A negative selection ensures T-cells do not recognize self-antigens. Here the remaining 10% of T-cells displaying high avidity for MHC bound peptides are eliminated by apoptosis. In many autoimmune diseases, this negative selection is defective, which leads to the production of autoreactive T cells.

Moving further down in the medulla, the DP T-cells recognize either MHC class I or MHC class II present on mTEC. When recognizing MHC-II, CD4 expression

is up-regulated whereas CD8 is down-regulated, giving birth to CD4+ T helper Lymphocytes. Conversely, when recognizing MHC-I, CD8 expression is up-regulated whereas CD4 is down-regulated, giving birth to CD8+ cytotoxic T-cells. In addition, agonist selection is a process enabling the deviation of intermediate avidity self-reactive CD4+ T-cells into thymic Tregs.

Thymic education is then completed, and T cells spread through our body, essentially in the peripheral lymph nodes where they encounter dendritic cells or other Antigen-Presenting Cells and continue their maturation and expansion.

B-cells’ primary development occurs in the bone marrow where hematopoietic stem cells successively differentiate into Pro B-cells, Pre-B cells, and finally immature (also called naïve) B-cells which carry a unique membrane-bound form of antibody constituting the antigen receptor or BCR. Immature B-cells move to the periphery e.g. spleen or lymph nodes for further maturation, activation, or differentiation into plasma cells.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



CELLULAR AND MOLECULAR MECHANISMS OF AI DISEASES

Central and peripheral tolerance

Central B-cell tolerance is controlled by multiple mechanisms at different B-cell differentiation stages, preventing the generation of auto-reactive B-cells and the subsequent autoantibody production involved in autoimmune diseases. In the bone marrow, Immature B cells encounter a wide variety of multivalent and monovalent self-antigens which are displayed by stromal cells, hematopoietic cells, and circulating molecules. If BCR from immature B cells do not, or only weakly, interact with multivalent self-antigens, they successfully leave the bone marrow and migrate to the peripheral lymphoid organs. Conversely, if the interaction between BCR and self-antigens is strong, further development of the potential auto-reactive B-cells is arrested. Three mechanisms are involved in acquiring central B-cell tolerance: receptor editing, clonal deletion, and anergy. Receptor editing is a process by which V and J genes encoding the BCR light chain undergo a new rearrangement. Receptor editing is the main process contributing to the elimination of autoreactive B cells. If receptor editing is not successful, then apoptosis of autoreactive B-cells occurs, which is known as clonal deletion. If BCR interact with monovalent soluble self-antigens, autoreactive immature B cells are functionally inactivated by a process called anergy. These anergic immature autoreactive B cells leave the bone marrow and migrate to the peripheral lymphoid organs, where they eventually undergo apoptosis. (54) (55) (56)

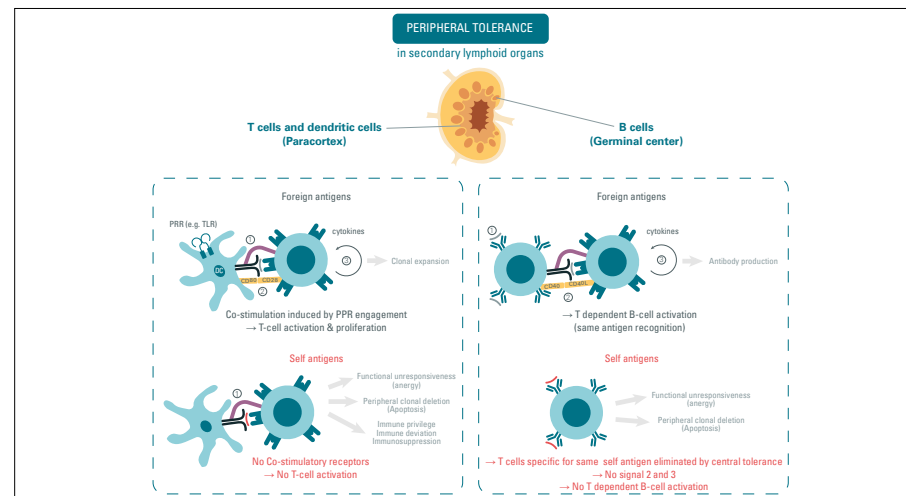
PERIPHERAL TOLERANCE

When central tolerance mechanisms are not efficient, self-reactive T and B cells escape and migrate to the peripheral lymphoid organs, where several peripheral tolerance mechanisms efficiently keep them in check. In fact, naïve T cells exhibit high levels of CCL21 receptor (CCR7) as well as receptors for adhesion molecules, and can only migrate to secondary lymphoid organs which specifically express CCL21 chemokine and adhesion molecules. Thus, naïve autoreactive T cells are excluded from peripheral tissues. In addition, some organs such as brain, eyes or testis are immune privileged areas where the immune cell numbers are intrinsically low (57).

Unlike DCs presenting foreign antigens, DCs presenting self-antigens express low levels of CD80 or CD86 co-stimulatory receptors which are necessary for complete T cell activation (58) (59). Thus, autoreactive T-cells recognizing MHC-II presenting

self-antigens cannot undergo activation but instead remain inactivated through anergy (60). Anergy, which refers to a long-term functional unresponsiveness state, can also result from the activation of inhibitory receptors such as CTLA-4, PD-1, and Lag3 present on the T-cell and which inhibit TCR signaling (61). Self-reactive T cells can undergo apoptosis through Activation Induced Cell Death (AICD) via Fas-Fas Ligand signaling and resulting in clonal deletion (62). Whereas clonal deletion and anergy prevent autoimmune responses, other mechanisms control or dampen autoimmune reactions. Among these peripheral tolerance mechanisms, Tregs are key players (see Regulatory T-cells in self-tolerance).

Autoreactive B cells which bind self-antigens in the periphery require co-stimulatory signals from helper T-cells (through CD40-CD40L and cytokines), which in such cases would be autoreactive T-cells. Thanks to the peripheral mechanisms cited above, autoreactive T-cells are eliminated. Thus, autoreactive B-cells which cannot receive activation signals from T-cells become anergic, and eventually die by apoptosis.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



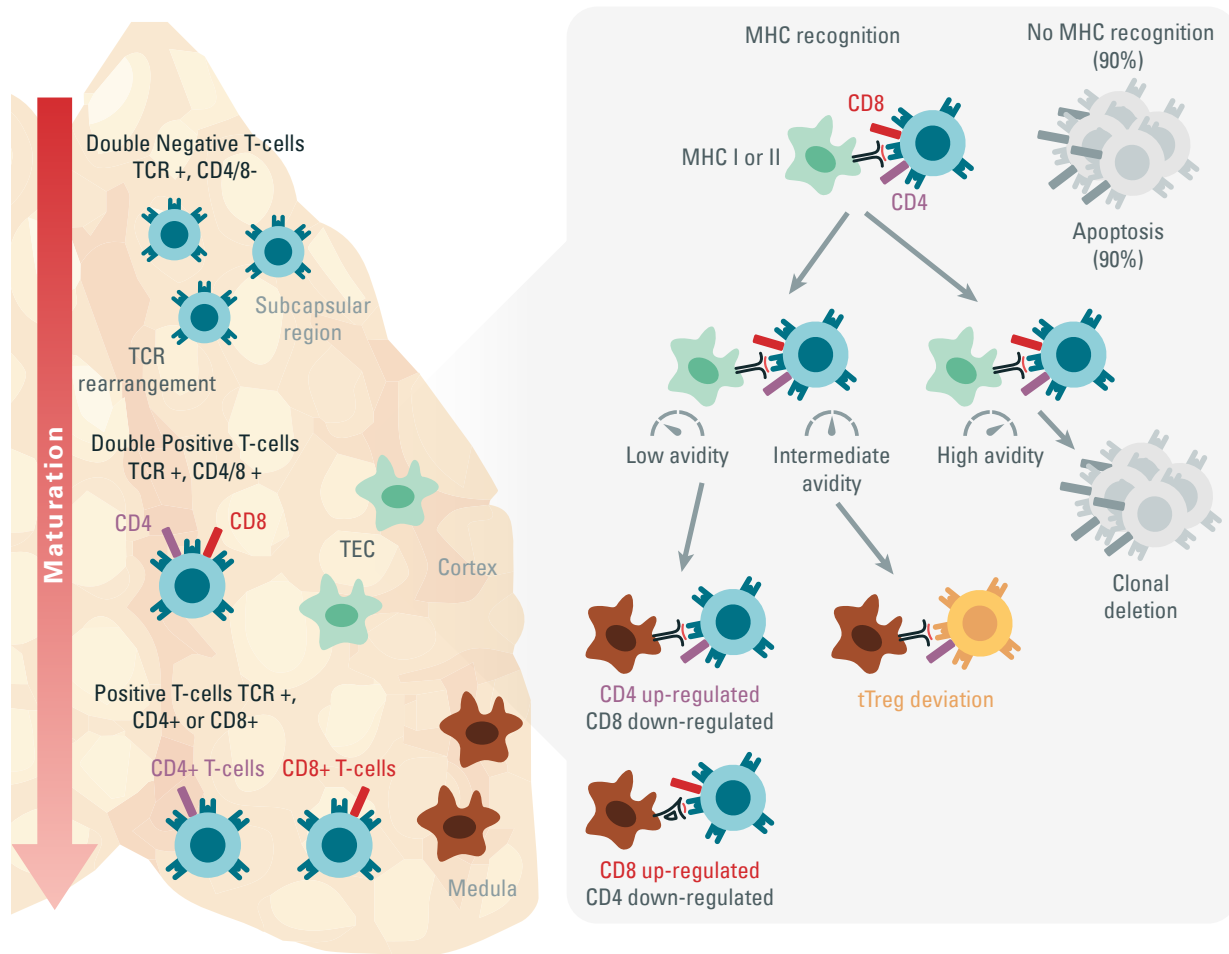
CENTRAL TOLERANCE

in primary lymphoid organs



Thymus: T cell selection & maturation

Normally, autoreactive T cells are eliminated early in their development but this process may be impaired in AI diseases



Bone marrow: B-cell maturation

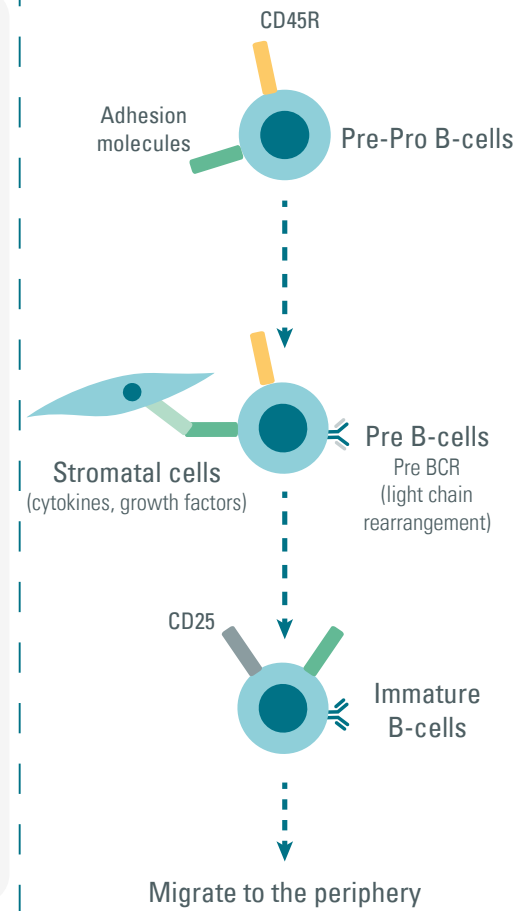


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



PERIPHERAL TOLERANCE

in secondary lymphoid organs

**T cells and dendritic cells
(Paracortex)**

**B cells
(Germinal center)**

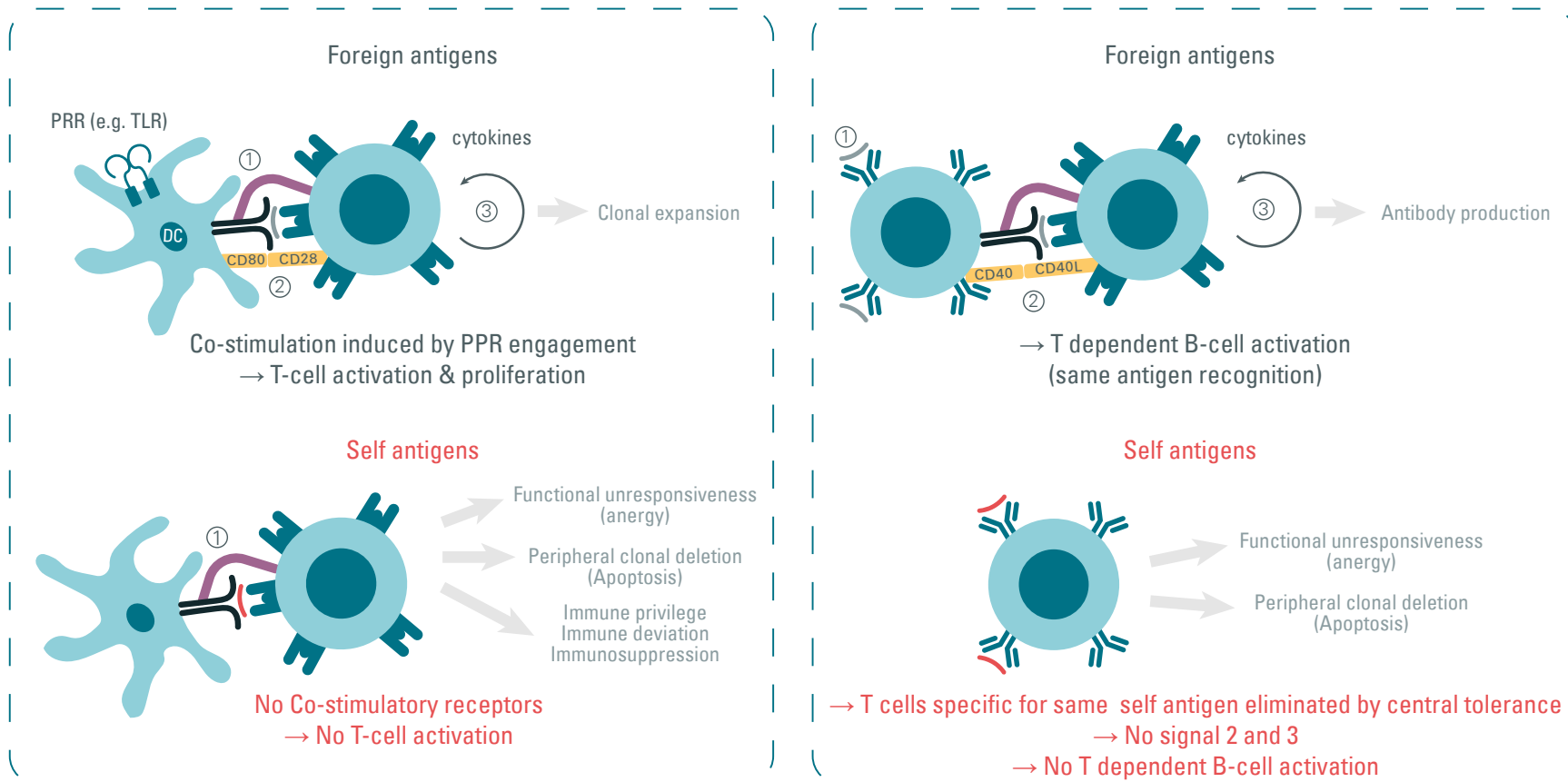
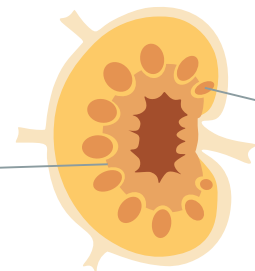


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

CELLULAR AND MOLECULAR MECHANISMS OF AI DISEASES

Central and peripheral tolerance

REGULATORY T-CELLS IN SELF TOLERANCE

If Thymic Tregs are involved in central T cell tolerance, peripheral Treg cells play essential roles in the immunosuppressive mechanisms leading to peripheral tolerance.

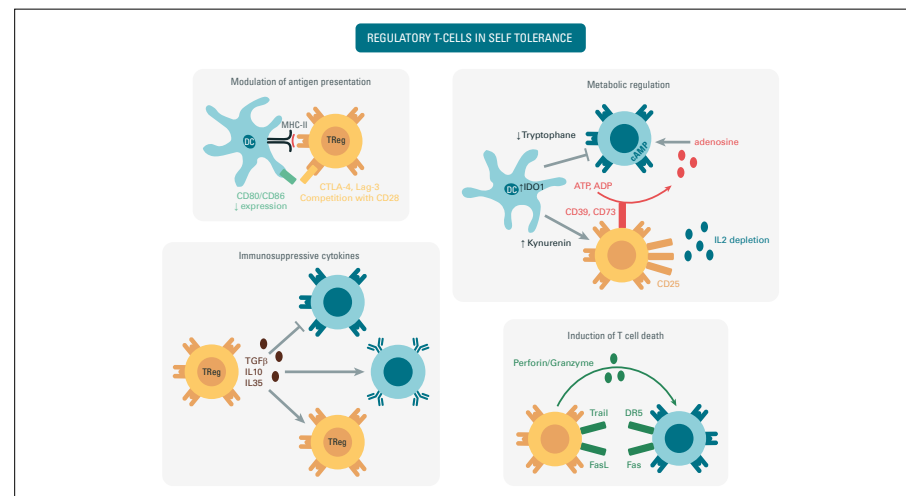
Direct and indirect immunosuppressive mechanisms exerted by Treg include:

- **Modulation of antigen presentation**, which is a key mechanism for Treg mediating suppression of effector T cells (63) (64). CTLA-4 is an immune checkpoint inhibitor constitutively expressed on the surface of Treg, which i) competes with CD28 expressed by effector T-cells for binding to CD80/86 ligands on DCs, and ii) down-regulates CD80/86 expression on DCs. Other immune checkpoint inhibitors like Lag-3 are also involved in this mechanism.
- **Metabolic regulation.** Tregs can induce the expression of Indoleamine 2,3-dioxygenase by DCs. Enzymatic activity of IDO catalyzes the production of kynurenine from Tryptophan, which promotes Treg differentiation through aryl hydrocarbon receptor (65) (66). In addition, Tregs express the CD39 and CD73 ectonucleotidases, which produce Adenosine from ATP or ADP. Adenosine binds the A2A receptor on effector T cells, which in turn results in their direct suppression due to increased levels of cAMP (67) (68).
- **Production of immunosuppressive cytokines.** Tregs can secrete TGFβ, IL10, or IL35, which inhibit effector T cells and DCs (69) (70) (71). In addition, Tregs express a high CD25 level which has high affinity for IL-2. Therefore, by consuming and depleting IL-2, Tregs prevent effector T cell activation (72).
- **Induction of effector T cell cytotoxicity.** Tregs can secrete Perforin, Granzyme A, and Granzyme B, which lead to effector T cells (73) (74). Apoptosis of Effector T cells induced by Tregs can be mediated by other molecules, such as TRAIL-DR5, FasL-Fas, Galectin-9-Tim3, and Gal-1 (75).

In addition to their role on effector T cells, Tregs are also involved in the suppression of autoantibody production by B cells (76).

Breakdown of central or peripheral tolerance mechanisms represents one cause of autoimmune disease onset.

It has been found that MHC gene variants are the most frequent genes associated with autoimmune diseases. Other genes related to immune defects, such as PTPN22 (protein phosphatase), CTLA4 (Checkpoint inhibitor), IL23R and TYK2, AIRE (transcription factor involved in central tolerance), FOXP3 (transcription factor involved in T-Reg lineage), IFIH1, DNASE1, TREX1, C1Q, or C4A (complement factors), have also been identified in certain autoimmune diseases.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



REGULATORY T-CELLS IN SELF TOLERANCE

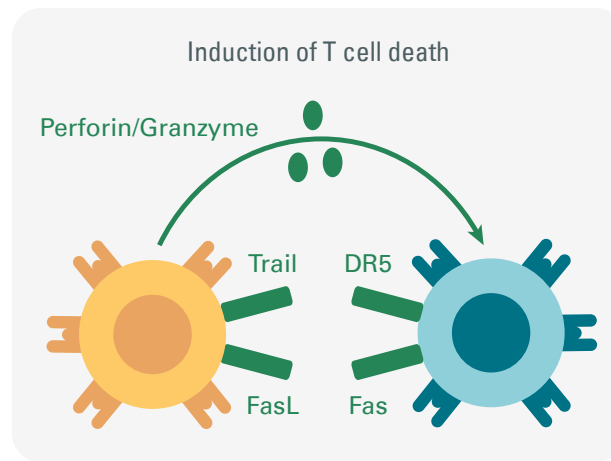
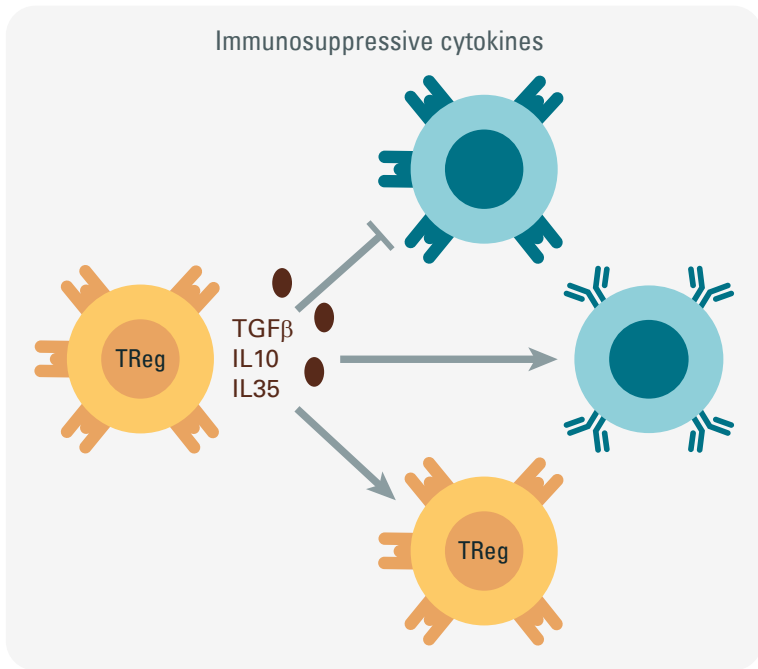
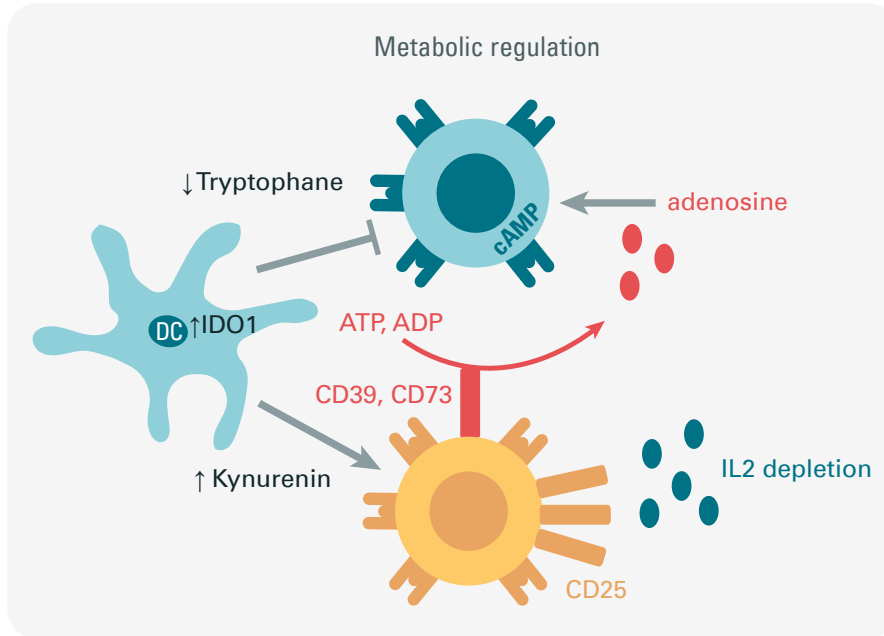
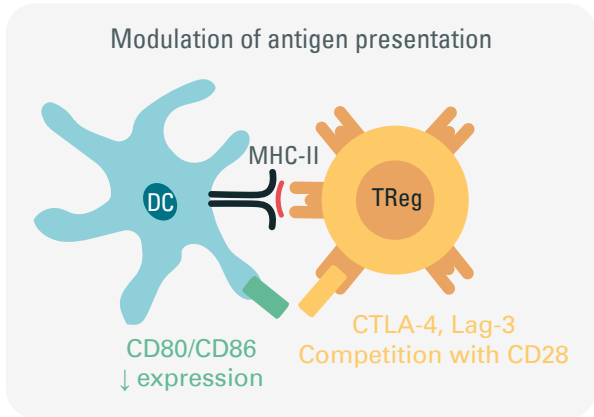


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

CELLULAR AND MOLECULAR MECHANISMS OF AI DISEASES

Innate immunity components

If a dysregulation of the adaptive immune response is involved in the development of autoimmune diseases, innate immune components which precede and prime the adaptive response also play a role in their pathophysiology.

Innate immune cells such as dendritic cells express a broad spectrum of sensors, called Pattern Recognition Receptors (PPR), that are either located at the plasma membrane or in the cytoplasm (77). PPRs recognize pathogen associated molecular patterns (PAMPs) and danger associated molecular patterns (DAMPs). Membrane receptors include Toll-like receptor 4, 1, or 2 (TLR), C-type lectin receptors (CLRs) and sialic acid binding Ig-like lectins (Siglecs). Endosomal sensors from the Toll-like receptor (TLR) family include TLR3 for double-strand RNA, TLR7 and TLR8 for single-strand RNA, and TLR9 for DNA (78) (79). Cytosolic sensors include Rig-1 like receptors (RLR) such as the helicases RIG-I for uncapped 5'-triphosphate RNA and MDA5 for long dsRNA (80); the cytosolic DNA Receptors (CDR) such as the double strand DNA cGAS enzyme or AIM2 (81), as well as cytosolic NOD-like Receptors (NLR) including NODs, NLRPs or NLRcs (82) (83). When activated, they all contribute to the transcriptional gene expression encoding proinflammatory cytokines such as TNF α , IL6, or IL1 β , and type 1 IFN. These cytokines play an essential role in bridging innate and adaptive immune responses. Activation of the TLRs, RLRs, or CDRs triggers signaling cascades which involve a variety of intracellular mediators such as Myd88, TRAF6, IRAKs, NFKB, or IRFs, and convey cytokine transcriptional programs.

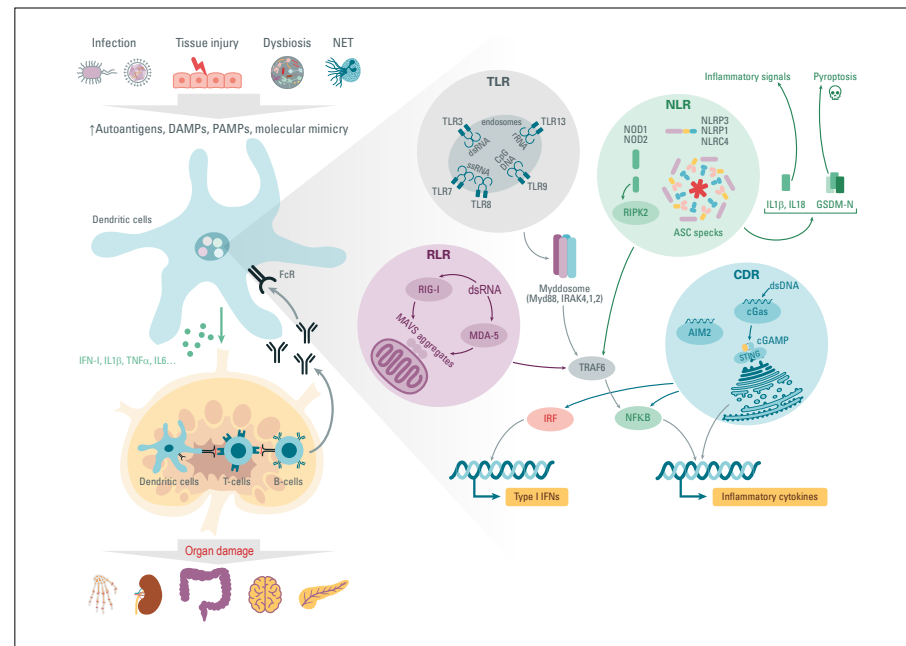
Inflammasome involves various NLR and contributes to the generation of the adaptive immune response through proteolytic maturation of inflammasome dependent cytokines (IL-1 β and IL-18). This enzymatic cleavage relies on ASC-speck formation and the subsequent Caspase-1 activation, and ultimately leads to an inflammatory programmed cell death called pyroptosis (84).

Numerous studies in animal and human models have revealed the crucial role of innate immune sensors in autoimmune pathogenesis. For example, lupus-mice models showed disease regression when IFNAR or TLR7 and 9 genes were deleted (85) (86). In human, the activation of MDA5 and RIG-1 has been reported in some SLE patients (87). Moreover, dysfunctions of DNase such as TREX1 contribute to various autoimmune diseases by the accumulation of DNA and subsequent PRR

activation (88). In addition, Neutrophil Extracellular Traps have been shown to induce TLR activation and type I IFN response (89). Inflammasome dysregulations have been associated with SLE and RA diseases (90). Furthermore, interaction between NETs and inflammasomes has been documented in SLE patients. SLE associated autoantibodies such as anti-dsDNA were shown to activate inflammasome in such patients (90).

Thus, innate immune components are now considered as major players in the innate immunity contribution to autoimmune pathogenesis.

For more information regarding general immunology and immune signaling pathways, please read our Immunology guide.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

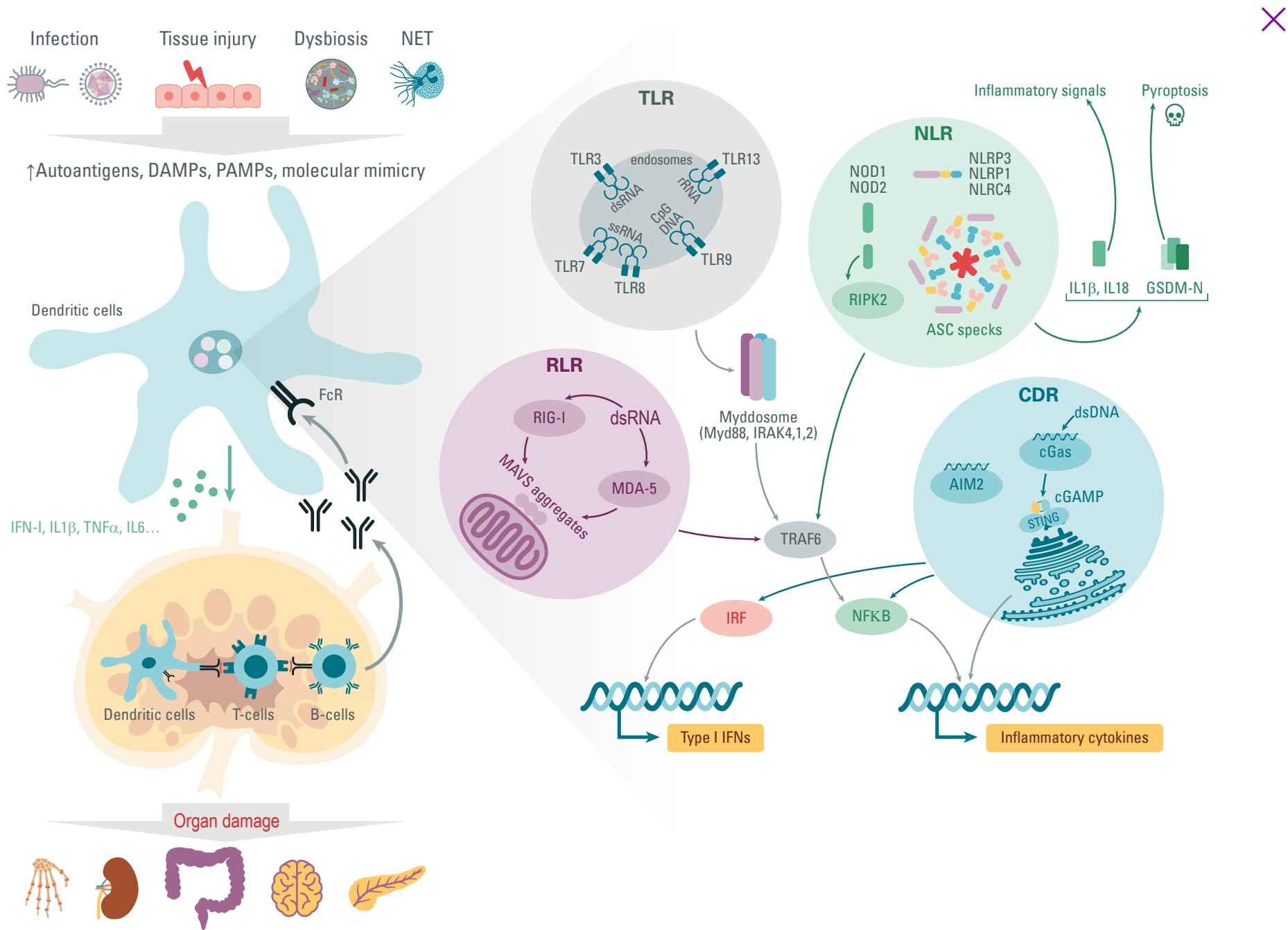


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

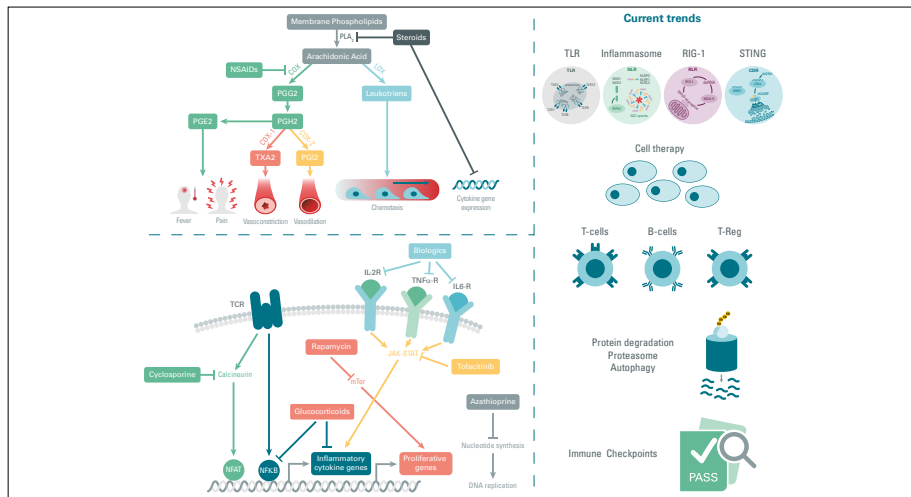
- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

THERAPEUTIC STRATEGIES IN AUTOIMMUNE DISEASE TREATMENT

Nonsteroidal anti-Inflammatory drugs(NSAIDs)

This class of molecules (e.g. ibuprofen) reduces inflammation by blocking the enzyme cyclooxygenases (COX-1 and COX-2) or COX-2 selectively (e.g. Celecoxib), and the production of thromboxane A2 and prostaglandins.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

THERAPEUTIC STRATEGIES IN AUTOIMMUNE DISEASE TREATMENT

Immunosuppressive drugs

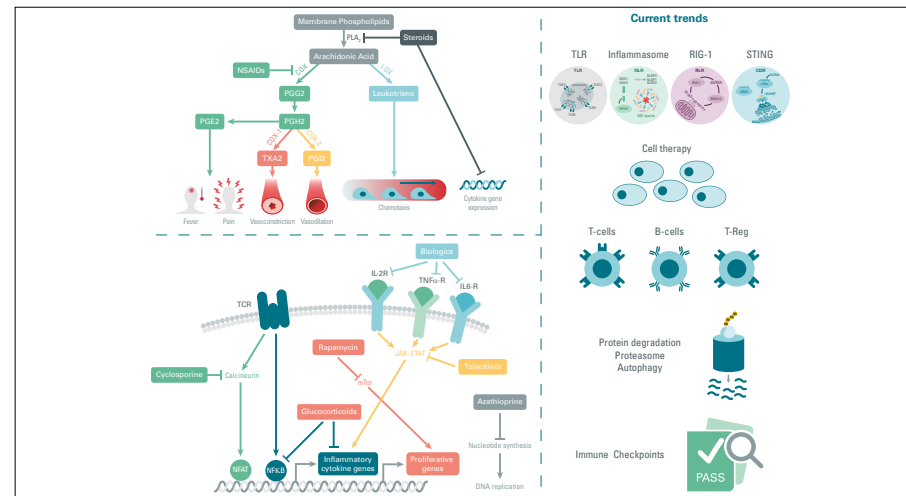
SMALL MOLECULES

- Glucocorticosteroids (e.g. cortisone) reduces inflammation by blocking Phospholipase A2 and the production of prostaglandins, leukotrienes and cytokines.
- Janus kinase inhibitors (e.g. tofacitinib) inhibit signal transduction from cytokine receptors.
- Calcineurin inhibitors (cyclosporine) inhibit the cytoplasmic phosphatase calcineurin, leading to reduced production of IL-2, IL-3, and interferon- γ .
- mTOR inhibitors (e.g. Rapamycin) inhibit intracellular FKBP-12 protein and reduce cytokine-induced T-cell proliferation.
- IMDH inhibitors (e.g. azathioprine) block nucleotide synthesis and lymphocyte proliferation.
- BTK inhibitors (e.g. Ibrutinib) inhibit BCR signaling and B-cell proliferation.

BIOLOGICS

- Anti-IL2 receptor antibodies (e.g. basiliximab, daclizumab) inhibit T cell proliferation by blocking the binding of IL-2.
- Anti-soluble TNF α antibodies (e.g. infliximab, Adalimumab) or fusion TNF receptor recombinant protein (etanercept) block TNF α pro inflammatory effects.
- Anti-IL6 receptor antibodies (e.g. tocilizumab) inhibit IL6 pro inflammatory effects.
- Anti-CD20 antibodies (e.g. Rituximab) induce B-cell apoptosis.

- Anti-CD49 (cell adhesion molecule α 4-integrin) antibodies (Natalizumab) block immune cell adhesion and their ability to pass through the intestinal barrier or blood brain barrier.
- Fusion CTLA-4 recombinant protein (Belatacept) inhibits T cell activation by blocking the B7 costimulatory molecule binding to CD28.
- Anti-VLA-4 antibodies (e.g. Natalizumab) block the VLA-4/VCAM-1 interaction and prevent immune cell infiltration.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

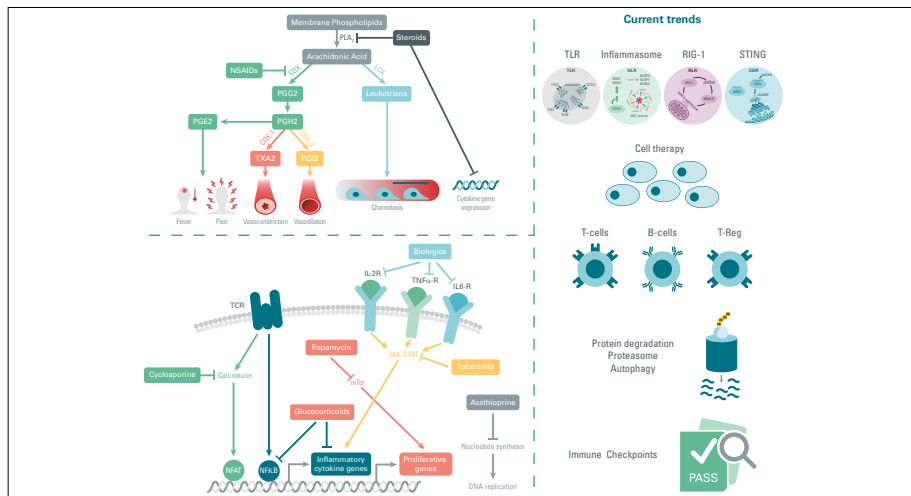
- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

THERAPEUTIC STRATEGIES IN AUTOIMMUNE DISEASE TREATMENT

Promising new therapeutic areas

Several new cellular and molecular targets have recently emerged as potential new areas for therapeutic intervention to treat auto-immune diseases. The targeting of innate immune components like inflammasome, TLR, or RLR is currently under scrutiny. Investigations aiming to restore immunosuppressive functions of T-Reg subsets or dampen B and T-cell activity may open up another area for future therapy, for example by targeting Immune checkpoint molecules. Intracellular pathways related to protein degradation (e.g. proteasome, autophagy) are being explored as well. Progress in cell therapies is also considered as a viable therapeutic axis in autoimmunity. Overall, improvements in autoimmune disease management may well result from the fast-growing understanding being achieved in the immunology field.



[Click to Enlarge](#)

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

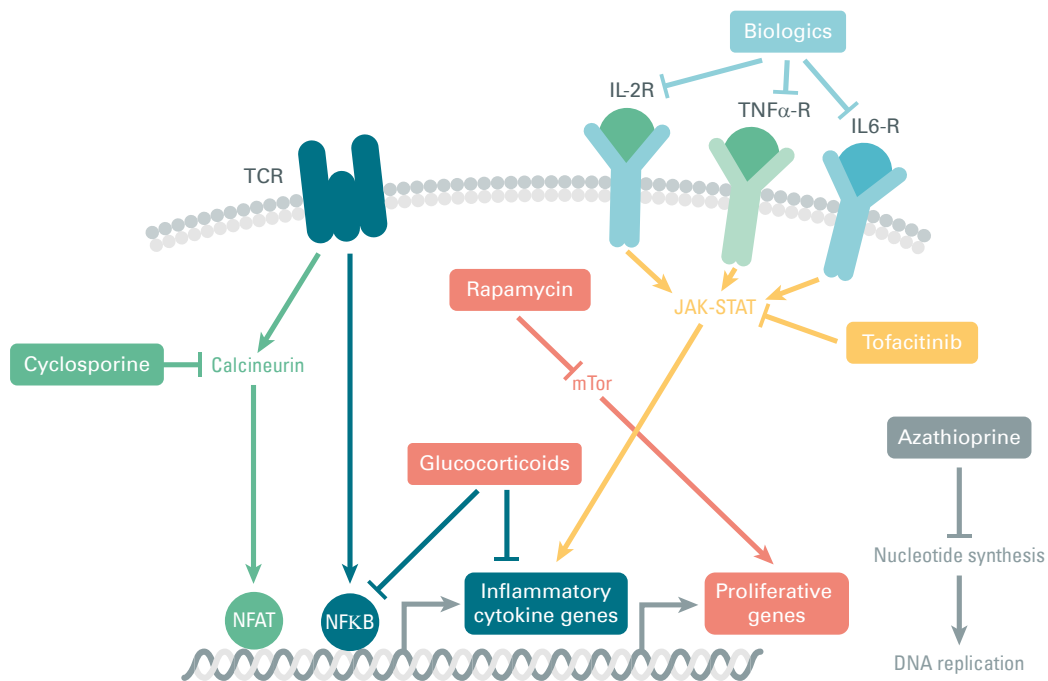
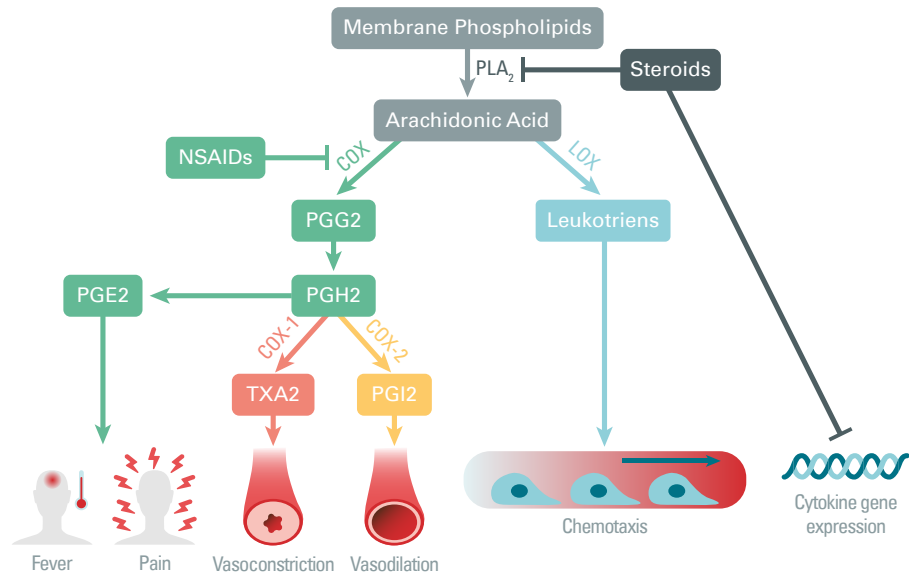
Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



Current trends

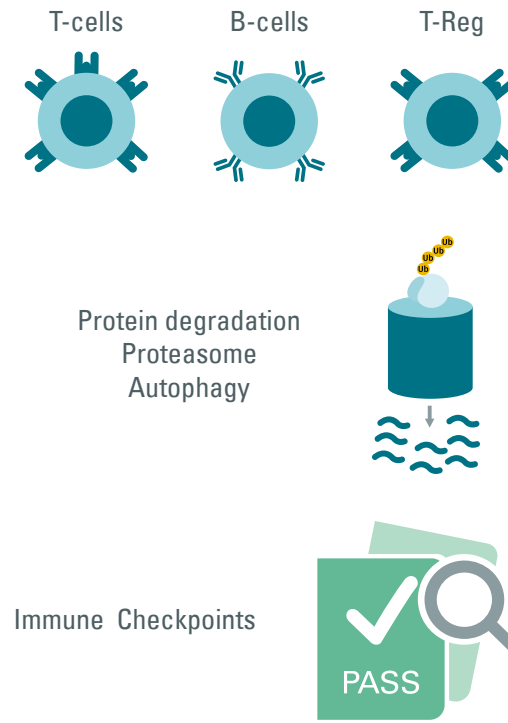
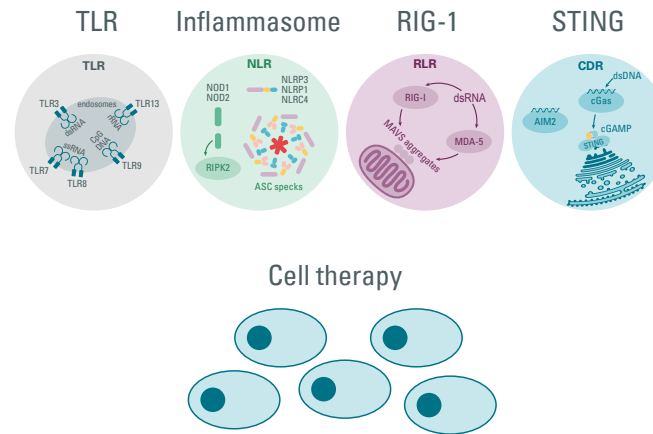


TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

BIBLIOGRAPHY

1. Incidence of rheumatoid arthritis in Sweden: a nationwide population-based assessment of incidence, its determinants, and treatment penetration. Eriksson JK, *et al.* s.l. : Arthritis Care Res, 2013, Vol. 65.
2. TK. Current tobacco smoking, formal education, and the risk of rheumatoid arthritis. Uhlig T, *et al.* s.l. : J Rheumatol., 1999, Vol. 26.
3. The contribution of HLA to rheumatoid arthritis. Deighton CM, *et al.* s.l. : Clin Genet., 1989, Vol. 36.
4. TRAF1-C5 as a risk locus for rheumatoid arthritis-a genome-wide study. Plenge RM, *et al.* s.l. : N Engl J Med. , 2007, Vol. 357.
5. STAT4 and the risk of rheumatoid arthritis and systemic lupus erythematosus. Remmers EF, *et al.* s.l. : N Engl J Med. , 2007, Vol. 357.
6. Epigenetics in the pathogenesis of RA. . Ospelt C, *et al.* s.l. : Semin Immunopathol., 2017, Vol. 39.
7. The epitopes targeted by the rheumatoid arthritis-associated antifilaggrin autoantibodies are posttranslationally generated on various sites of (pro)filaggrin by deimination of arginine residues. Girbal-Neuhauser E, *et al.* s.l. : J Immunol., 1999, Vol. 162.
8. The American College of Rheumatology / European League Against Rheumatism Classification Criteria for Rheumatoid Arthritis. Aletaha D, *et al.* s.l. : Ann Rheum Dis., 2010, Vol. 69.
9. Cytokines in rheumatoid arthritis - shaping the immunological landscape. McInnes, I., *et al.* s.l. : Nat Rev Rheumatol., 2016, Vol. 12.
10. The worldwide incidence and prevalence of systemic lupus erythematosus: a systematic review of epidemiological studies. Rees F, *et al.* s.l. : Rheumatology., 2017, Vol. 56.
11. Complement deficiencies and systemic lupus erythematosus. Truedsson, L. *et al.* s.l. : Autoimmunity, 2007, Vol. 40.
12. Risk alleles for systemic lupus erythematosus in a large case-control collection and associations with clinical subphenotypes. Taylor, K. E. *et al.* s.l. : PLoS Genet., 2011, Vol. 7.
13. Mutations in the gene encoding the 3'-5' DNA exonuclease TREX1 are associated with systemic lupus erythematosus. Lee-Kirsch MA, *et al.* s.l. : Nat Genet., 2007, Vol. 39.
14. Estrogen and autoimmune disease. SE., Walker. s.l. : Clin Rev Allergy Immunol. , 2011, Vol. 40.
15. Systemic lupus erythematosus. Kaul, A., *et al.* s.l. : Nat Rev Dis Primers, 2016.
16. Neutrophil extracellular traps kill bacteria. Brinkmann V, *et al.* s.l. : Science, 2004, Vol. 303.
17. Nucleic acids of mammalian origin can act as endogenous ligands for Toll-like receptors and may promote systemic lupus erythematosus. Barrat, F. J. *et al.* 2005 : J. Exp. Med., Vol. 202.
18. Dysregulation of antiviral helicase pathways in systemic lupus erythematosus. Oliveira, L. *et al.* s.l. : Front. Genet., 2014, Vol. 5.
19. Autoantigen microarray for high throughput autoantibody profiling in systemic lupus erythematosus. Zhu, H. *et al.* s.l. : Genomics Proteomics Bioinformatics, 2015, Vol. 13.
20. T cells as therapeutic targets in SLE. Crispin, J. C., *et al.* s.l. : Nat. Rev. Rheumatol., 2010, Vol. 6.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

BIBLIOGRAPHY

21. Expansion of circulating T cells resembling follicular helper T cells is a fixed phenotype that identifies a subset of severe systemic lupus erythematosus. Simpson, N. *et al.* Arthritis Rheum. 2010, Vol. 62.
22. Autoantibody explosion in systemic lupus erythematosus: erythematosus: more than 100 different antibodies found in SLE patients. Sherer Y, *et al.* s.l. : Semin Arthritis Rheum., 2004, Vol. 34.
23. Long-lived autoreactive plasma cells drive persistent autoimmune inflammation. Hiepe, F. *et al.* s.l. : Nat. Rev. Rheumatol., 2011, Vol. 7.
24. Association analyses identify 38 susceptibility loci for inflammatory bowel disease and highlight shared genetic risk across populations. Liu, J., *et al.* s.l. : Nat Genet., 2015, Vol. 47.
25. A genome-wide association scan of nonsynonymous SNPs identifies a susceptibility variant for Crohn's Disease in ATG16L1. Hampe J, *et al.* s.l. : Nat Genet., 2007, Vol. 39.
26. Epidemiology of inflammatory bowel disease. Loftus Jr EV, Sandborn WJ. s.l. : Gastroenterol Clin North Am. , 2002, Vol. 31.
27. The genetic background of inflammatory bowel disease: from correlation to causality. Uniken Venema WT, *et al.* s.l. : J Pathol., 2017, Vol. 241.
28. The microbiota in adaptive immune homeostasis and disease. Honda, K., *et al.* s.l. : Nature, 2016, Vol. 535.
29. The IL-23-IL-17 immune axis: from mechanisms to therapeutic testing. 45. Gaffen SL, Jain R, Garg AV, Cua DJ. s.l. : Nat Rev Immunol. , 2014, Vol. 14.
30. IL-1 β mediates chronic intestinal inflammation by promoting the accumulation of IL-17A secreting innate lymphoid cells and CD4+ Th17 cells. Coccia, M., *et al.* s.l. : J Exp Med, 2012, Vol. 209.
31. Neutralisation of the interleukin-33/ST2 pathway ameliorates experimental colitis through enhancement of mucosal healing in mice. Sedhom MAK, *et al.* s.l. : Gut, 2013, Vol. 62.
32. ASCA, ANCA, ALCA and many more: are they useful in the diagnosis of inflammatory bowel disease? . Zhou G, *et al.* s.l. : Dig Dis. , 2016;
33. Pleiotropic functions of TNF- α in the regulation of the intestinal epithelial response to inflammation. Leppkes M, *et al.* s.l. : Int Immunol., 2014, Vol. 26.
34. Multiple sclerosis. Noseworthy JH, *et al.* s.l. : N Engl J Med., 2000, Vol. 343.
35. Interactions between genetic, lifestyle and environmental risk factors for multiple sclerosis. . Olsson T, *et al.* s.l. : Nat Rev Neurol., 2017, Vol. 13.
36. HLA-A confers an HLA-DRB1 independent influence on the risk of multiple sclerosis. . Brynedal B, *et al.* s.l. : PLoS One., 2007.
37. Epistasis among HLA-DRB1, HLA-DQA1, and HLA-DQB1 loci determines multiple sclerosis susceptibility. Lincoln MR, *et al.* s.l. : Proc Natl Acad Sci U.S.A., 2009, Vol. 106.
38. Identification of Epstein-Barr virus proteins as putative targets of the immune response in multiple sclerosis. Cepok S, *et al.* s.l. : J Clin Invest, 2005, Vol. 115.
39. Environmental and genetic risk factors for MS: an integrated review. Waubant, E. *et al.* s.l. : Annals of Clinical and Translational Neurology, 2019, Vol. 6.
40. molecular mimicry in multiple sclerosis. Libbey J, *et al.* s.l. : International review of neurobiology, 2007, Vol. 79.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography



BIBLIOGRAPHY

41. Current multiple sclerosis treatments have improved our understanding of MS autoimmune pathogenesis. *al., Martin R et. s.l. : Eur. J. Immunol., 2016, Vol. 46.*
42. Immunoblot detection of oligoclonal anti-myelin basic protein IgG antibodies in cerebrospinal fluid in multiple sclerosis. *Cruz M, et al. s.l. : Neurology., 1987, Vol. 37.*
43. A study of immunoglobulin G in the cerebrospinal fluid of 1007 patients with suspected neurological disease using isoelectric focusing and the Log IgG-Index
A comparison and diagnostic applications. *Mclean BN, et al. s.l. : Brain. , 1990, Vol. 113.*
44. Conformational epitopes of myelin oligodendrocyte glycoprotein are targets of potentially pathogenic antibody responses in multiple sclerosis. *Menge T, et al. s.l. : J Neuroinflamm., 2011, Vol. 8.*
45. Staging presymptomatic type 1 diabetes: a scientific statement of JDRF, the Endocrine Society, and the American Diabetes Association. *Insel RA, et al. s.l. : Diabetes Care., 2015, Vol. 38.*
46. Genetic risk factors for type 1 diabetes. *Pociot F, Lernmark A. s.l. : Lancet., 2016, Vol. 387.*
47. Type 1 diabetes mellitus. *Katsarou A, et al. s.l. : Nat Rev Dis Primers., 2017, Vol. 3.*
48. Patterns of β -cell autoantibody appearance and genetic associations during the first years of life. *Ilonen, J. et al. s.l. : Diabetes , 2013, Vol. 62.*
49. V(D)J recombination and the evolution of the adaptive immune system. *Market E, Papavasiliou FN. s.l. : PLoS Biol., 2003.*
50. Cell death and thymic tolerance. *Daley SR, et al. s.l. : Immunol Rev., 2017, Vol. 277.*
51. Positive selection of thymocytes. . *Jameson SC, Hogquist KA, Bevan MJ. s.l. : Annu Rev Immunol., 1995, Vol. 13.*
52. Negative selection – clearing out the bad apples from the T-cell repertoire. . *E., Palmer. s.l. : Nat Rev Immunol. , 2003, Vol. 3.*
53. Positive and negative selection of the T cell repertoire: what thymocytes see (and don't see). *Klein L, Kyewski B, Allen PM, Hogquist KA. s.l. : Nat Rev Immunol., 2014, Vol. 14.*
54. Receptor editing is the main mechanism of B cell tolerance toward membrane antigens. *Halverson R, Torres RM, Pelanda R. s.l. : Nat Immunol., 2004, Vol. 5.*
55. Central B cell tolerance: where selection begins. *Pelanda R, Torres RM. s.l. : Cold Spring Harb Perspect Biol., 2012, Vol. 4.*
56. The scope of receptor editing and its association with autoimmunity. . *Verkoczy LK, Martensson AS, Nemazee D. s.l. : Curr Opin Immunol., 2004, Vol. 16.*
57. What is immune privilege (not)? *Galea I, Bechmann I, Perry VH. s.l. : Trends Immunol., 2007, Vol. 28.*
58. Antigen-specific inhibition of effector T cell function in humans after injection of immature dendritic cells. *Dhodapkar MV, et al. 193, s.l. : J Exp Med., 2001, Vol. 2001.*
59. Recent thymic emigrants are tolerized in the absence of inflammation. . *Friesen TJ, et al. s.l. : J Exp Med., 2016, Vol. 213.*
60. T-cell unresponsiveness *in vivo* and *in vitro*: fine specificity of induction and molecular characterization of the unresponsive state. *Jenkins MK, et al. s.l. : Immunol Rev., 1987, Vol. 95.*

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

BIBLIOGRAPHY

61. NFAT, immunity and cancer: a transcription factor comes of age. Müller MR, Rao A. s.l. : Nat Rev Immunol., 2010, Vol. 10.
62. Fas and the art of lymphocyte maintenance. MJ., Lenardo. s.l. : J Exp Med., 1996, Vol. 183.
63. CTLA-4 control over Foxp3+ regulatory T cell function. Wing K, Onishi Y, Prieto-martin P, *et al.* s.l. : Science., 2008, Vol. 322.
64. Regulatory T cells inhibit dendritic cells by lymphocyte activation gene-3 engagement of MHC class II. Liang B, Workman C, Lee J, *et al.* s.l. : J Immunol., 2008, Vol. 180.
65. Modulation of tryptophan catabolism by regulatory T cells. . Fallarino F, Grohmann U, Hwang KW, *et al.* s.l. : Nat Immunol., 2003, Vol. 4.
66. An interaction between kynurenine and the aryl hydrocarbon receptor can generate regulatory T cells. Mezrich JD, *et al.* s.l. : J Immunol., 2010, Vol. 185.
67. T regulatory and primed uncommitted CD4 T cells express CD73, which suppresses effector CD4 T cells by converting 5'-adenosine monophosphate to adenosine. Kobie JJ, *et al.* s.l. : J Immunol., 2006, Vol. 177.
68. Adenosine generation catalyzed by CD39 and CD73 expressed on regulatory T cells mediates immune suppression. Deaglio S, *et al.* s.l. : J Exp Med., 2007, Vol. 204.
69. Cell contact-dependent immunosuppression by CD4(+)CD25(+) regulatory T cells is mediated by cell surface-bound transforming growth factor beta. Nakamura K, *et al.* s.l. : J Exp Med., 2001, Vol. 194.
70. An essential role for interleukin 10 in the function of regulatory T cells that inhibit intestinal inflammation. Asseman C, *et al.* s.l. : J Exp Med., 1999, Vol. 190.
71. IL-35-mediated induction of a potent regulatory T cell population. Collison LW, *et al.* s.l. : Nat Immunol., 2010, Vol. 11.
72. Interleukin-2 is essential for CD4+CD25+ regulatory T cell function. de la Rosa M, *et al.* s.l. : Eur J Immunol., 2004, Vol. 34.
73. Differential expression of granzymes A and B in human cytotoxic lymphocyte subsets and T regulatory cells. Grossman WJ, *et al.* s.l. : Blood., 2004, Vol. 104.
74. Granzyme B and perforin are important for regulatory T cell-mediated suppression of tumor clearance. Cao X, *et al.* s.l. : Immunity. , 2007, Vol. 27.
75. Tim-3-Galectin-9 pathway involves the suppression induced by CD4+CD25+ regulatory T cells. Wang F, *et al.* s.l. : Immunobiology, 2009, Vol. 214.
76. CD4+ CD25+ regulatory T cells inhibit the maturation but not the initiation of an autoantibody response. Fields ML, *et al.* s.l. : J Immunol., 2005, Vol. 175.
77. Pattern recognition receptors and inflammation. Takeuchi O, Akira S. s.l. : Cell., 2010, Vol. 140.
78. The history of Toll-like receptors – redefining innate immunity. O'Neill LA, Golenbock D, Bowie AG. s.l. : Nat Rev Immunol., 2013, Vol. 13.
79. The role of pattern-recognition receptors in innate immunity: update on Toll-like receptors. . Kawai T, Akira S. s.l. : Nat Immunol., 2010, Vol. 11.
80. Viral RNA detection by RIG-I-like receptors. Yoneyama M, *et al.* s.l. : Curr Opin Immunol., 2015, Vol. 32.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
- Systemic lupus erythematosus
- Inflammatory bowel disease
- Multiple sclerosis
- Type 1 diabetes
- Common antigens and autoantibodies associated with AI

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
- Innate immunity components

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
- Immunosuppressive drugs
- Promising new therapeutic areas

Bibliography

BIBLIOGRAPHY

81. Intracellular nucleic acid detection in autoimmunity. Crowl JT, *et al.* s.l. : Annu Rev Immunol., 2017, Vol. 35.
82. NOD1 and NOD2: signaling, host defense, and inflammatory disease. Caruso R, *et al.* s.l. : Immunity., 2014, Vol. 41.
83. Differential requirement for Caspase-1 autoproteolysis in pathogen-induced cell death and cytokine processing. Broz P, *et al.* s.l. : Cell Host Microbe., 2010, Vol. 8.
84. Converging roles of caspases in inflammasome activation, cell death and innate immunity. . Man SM, Kanneganti TD. s.l. : Nat Rev Immunol., 2016, Vol. 16.
85. Type-I interferon receptor deficiency reduces lupus-like disease in NZB mice. Santiago-Raber, M.L. *et al.* s.l. : J. Exp. Med., 2003, Vol. 197.
86. Toll-like receptor 7 and TLR9 dictate autoantibody specificity and have opposing inflammatory and regulatory roles in a murine model of lupus. Christensen, S.R. *et al.* s.l. : Immunity, 2006, Vol. 25.
87. Reactive oxygen species induce virus-independent MAVS oligomerization in systemic lupus erythematosus. Buskiewicz, I.A. *et al.* s.l. : Sci. Signal., 2016.
88. Theofilopoulos, A.N., *et al.* Intracellular nucleic acid sensors and autoimmunity. s.l. : J. Interferon Cytokine Res., 2011, Vol. 31.
89. Neutrophils activate plasmacytoid dendritic cells by releasing self-DNA-peptide complexes in systemic lupus erythematosus. Lande, R. *et al.* s.l. : Sci. Transl. Med., 2011.
90. Inflammasomes and autoimmune and rheumatic diseases: A comprehensive review. J.I. Shin, *et al.* s.l. : Journal of Autoimmunity 103, 2019, Vol. 103.
91. NFAT,immunity and cancer: a transcription factor comes of age. Müller MR, Rao A. s.l. : Nat Rev Immunol., 2010, Vol. 10.

TABLE OF CONTENTS

Purpose and scope

Autoimmune diseases at a glance

Pathophysiology of autoimmune diseases

- Rheumatoid arthritis
 - Systemic lupus erythematosus
 - Inflammatory bowel disease
 - Multiple sclerosis
 - Type 1 diabetes
 - Common antigens and autoantibodies associated with AI
-

Cellular and molecular mechanisms of AI diseases

- Central and peripheral tolerance
 - Innate immunity components
-

Therapeutic strategies in autoimmune disease treatment

- Nonsteroidal anti-inflammatory drugs (NSAIDs)
 - Immunosuppressive drugs
 - Promising new therapeutic areas
-

Bibliography



www.revivity.com

revivity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
(800) 762-4000 | www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2023, Revvity, Inc. All rights reserved.

203053