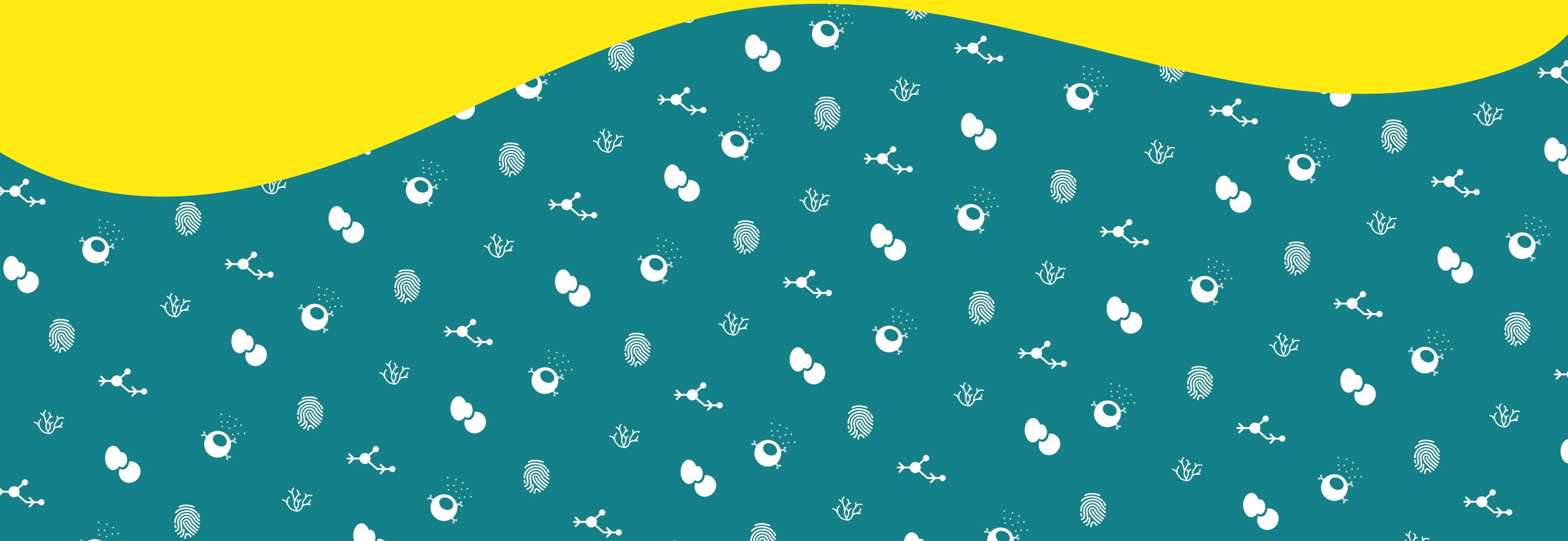


revvity

HTRF solutions

Guide to major applications



AUTOIMMUNITY

HTRF solutions for autoimmunity

Revvity's kits and services offer simple, rapid, cost-effective solutions for auto-immunity research.

Referenced in a growing number of publications aiming at understanding the pathophysiology of Rheumatoid Arthritis, Lupus, Multiple Sclerosis and Inflammatory Bowel Disease.

Comprehensive portfolio of Total and Phospho-protein, cytokine, chemokine and kinase assays, covering all major pro-inflammatory pathways, including **NF-κB**, **ZAP-70**, **STING**, **NLRP3**, and **JAK-STAT**.

	Specific cytokines of interest
Lupus	<i>IFN alpha, beta, gamma</i>
Psoriasis	<i>TNF alpha, IL8, IL17</i>
Multiple Sclerosis	<i>TNF alpha</i>
Inflammatory Bowel Disease	<i>TNF alpha, IL6, IL23</i>
Autoimmune arthritis	<i>TNF alpha, IL6, IL17, IL23, GM-CSF</i>

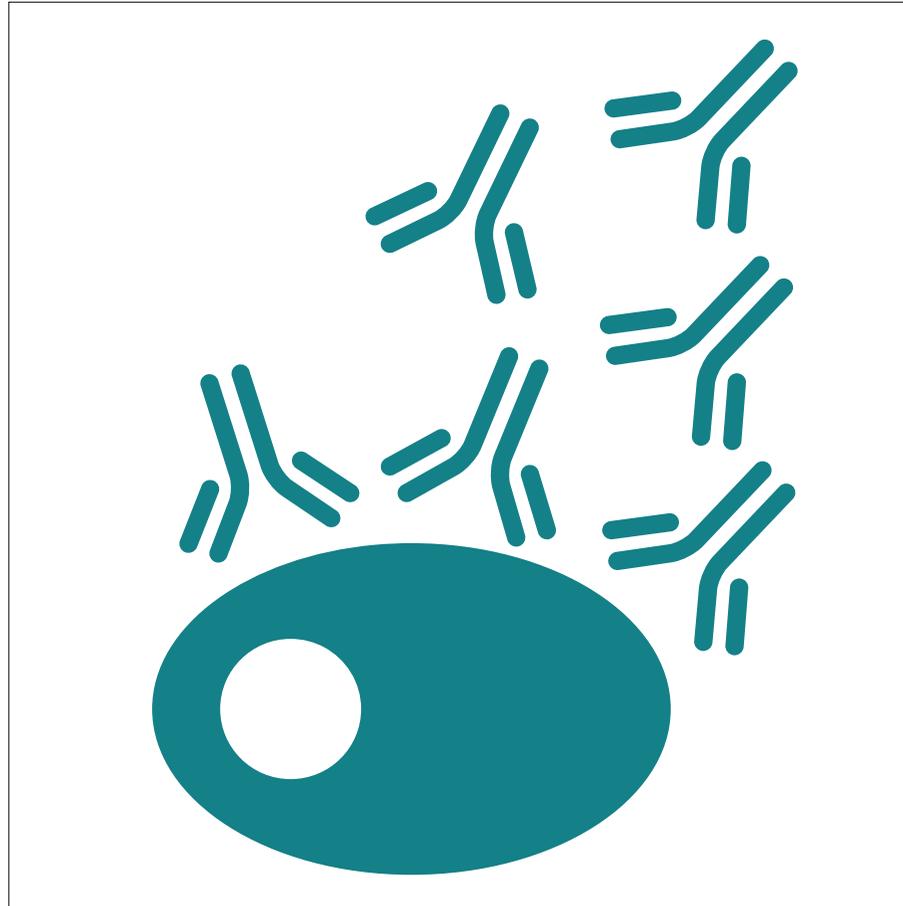


TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)

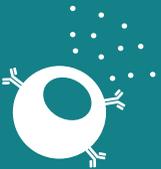
DISEASES	Lupus Multiple Sclerosis	Lupus Psoriasis Multiple Sclerosis Inflammatory Bowel Disease Rheumatoid arthritis	Lupus Autoimmune arthritis	Lupus	Lupus Psoriasis Multiple Sclerosis Inflammatory Bowel Disease Rheumatoid arthritis
	STING	JAK-STAT	BTK	ZAP70	NF-kB
Cytokines 	<i>IFN beta</i> <i>IFN alpha</i> <i>IFN gamma</i>	<i>IFN beta</i> <i>IFN alpha</i>	<i>IFN gamma</i> <i>TNF alpha</i> <i>IL2</i> <i>IL4</i> <i>IL5</i> <i>IL6</i> <i>IL10</i> <i>IL12</i> <i>IL17</i>	<i>TNF alpha</i> <i>IFN gamma</i> <i>IL2</i> <i>IL4</i> <i>IL5</i> <i>IL6</i> <i>IL10</i>	<i>IFN alpha</i> <i>IFN beta</i> <i>IFN gamma</i> <i>IL6</i> <i>CCL2</i> <i>CCL3</i> <i>CCL4</i> <i>CCL5</i> <i>CCL20</i> <i>CXCL1</i> <i>CXCL10</i>
Phospho-proteins 	<i>Alpha Tubulin Total & Phospho-STING (Ser366)</i> <i>Phospho-IRF3 (Ser386) Total & Phospho-TBK1 (Ser172)</i> <i>Total & Phospho-NF-kB (Ser536)</i> <i>Total & Phospho-STAT1 (Tyr701)</i> <i>Total & Phospho-STAT3 (Tyr705)</i> <i>Total & Phospho-STAT5 (Tyr694)</i> <i>Total IRAK3</i> <i>Total IRAK4</i>	<i>Alpha Tubulin Total & Phospho-STAT1 (Tyr701)</i> <i>Total & Phospho-STAT3 (Tyr705)</i> <i>Total & Phospho-STAT5 (Tyr694)</i> <i>Total & Phospho-AKT1 (Ser473)</i> <i>Total & Phospho-AKT2 (Ser473)</i> <i>Total & Phospho-AKT3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Thr308)</i> <i>Total AKT1/2/3 (Ser2448)</i> <i>Phospho-mTOR (Ser2448)</i> <i>Total & Phospho-ERK (Thr202/Tyr204)</i>	<i>Alpha Tubulin Total & Phospho-BTK (Tyr223)</i> <i>Total & Phospho-AKT1 (Ser473)</i> <i>Total & Phospho-AKT2 (Ser473)</i> <i>Total & Phospho-AKT3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Thr308)</i> <i>Total AKT1/2/3 (Ser2448)</i> <i>Phospho-mTOR (Ser2448)</i> <i>Total & Phospho-ERK (Thr202/Tyr204)</i> <i>Phospho-JNK (Thr183/Tyr185)Total & Phospho-NF-kB (Ser536)</i> <i>Total & Phospho-GSK3 (Ser9)</i> <i>Phospho-BAD (ser112)</i> <i>Phospho-P38 (thr180/Tyr182)</i>	<i>Alpha Tubulin Total & Phospho-ZAP70 (Tyr319)</i> <i>Total & Phospho-SLP-76 (Ser376)</i> <i>Total & Phospho-CREB (Ser133)</i> <i>Total & Phospho-ERK (Thr202/Tyr204)</i> <i>Phospho-JNK (Thr183/Tyr185)Total & Phospho-AKT1 (Ser473)</i> <i>Total & Phospho-AKT2 (Ser473)</i> <i>Total & Phospho-AKT3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Thr308)</i> <i>Total AKT1/2/3 (Ser2448)</i> <i>Phospho-mTOR (Ser2448)</i> <i>Total & Phospho-SHP1 (Tyr564)</i> <i>Total & Phospho-SHP2 (Tyr542)</i>	<i>Alpha Tubulin Total & Phospho-NF-kB (Ser536)</i> <i>Total & Phospho-IKKbeta (Ser177/181)</i> <i>Total IRAK3</i> <i>Total IRAK4</i>

TABLE OF CONTENTS

- [HTRF solutions for autoimmunity](#)
- [HTRF solutions for cardiology](#)
- [HTRF solutions for NAFLD and fibrosis](#)
- [HTRF solutions for neuroscience](#)
- [HTRF solutions for CAR-T research](#)
- [HTRF solutions for TLR research](#)
- [HTRF solutions for virology](#)



DISEASES	Lupus Multiple Sclerosis	Lupus Psoriasis Multiple Sclerosis Inflammatory Bowel Disease Rheumatoid arthritis	Lupus Autoimmune arthritis	Lupus	Lupus Psoriasis Multiple Sclerosis Inflammatory Bowel Disease Rheumatoid arthritis
	STING	JAK-STAT	BTK	ZAP70	NF-kB
 Kinases (Biochemical)		JAK1 (TK) JAK2 (TK) JAK3 (TK)	LYN (TK) SYK (TK)	LCK (TK)	
 PPI	PD1/PDL1 CTLA4/B7-1 CTLA4/B7-2 LAG3/MHCII Human H232 STING binding kit Human AQ STING binding kit Mouse STING binding kit	CD47/SIRPa		PD1/PDL1 CTLA4/B7-1 CTLA4/B7-2	

TABLE OF CONTENTS

HTRF solutions for autoimmunity

HTRF solutions for cardiology

HTRF solutions for NAFLD and fibrosis

HTRF solutions for neuroscience

HTRF solutions for CAR-T research

HTRF solutions for TLR research

HTRF solutions for virology

CARDIOLOGY

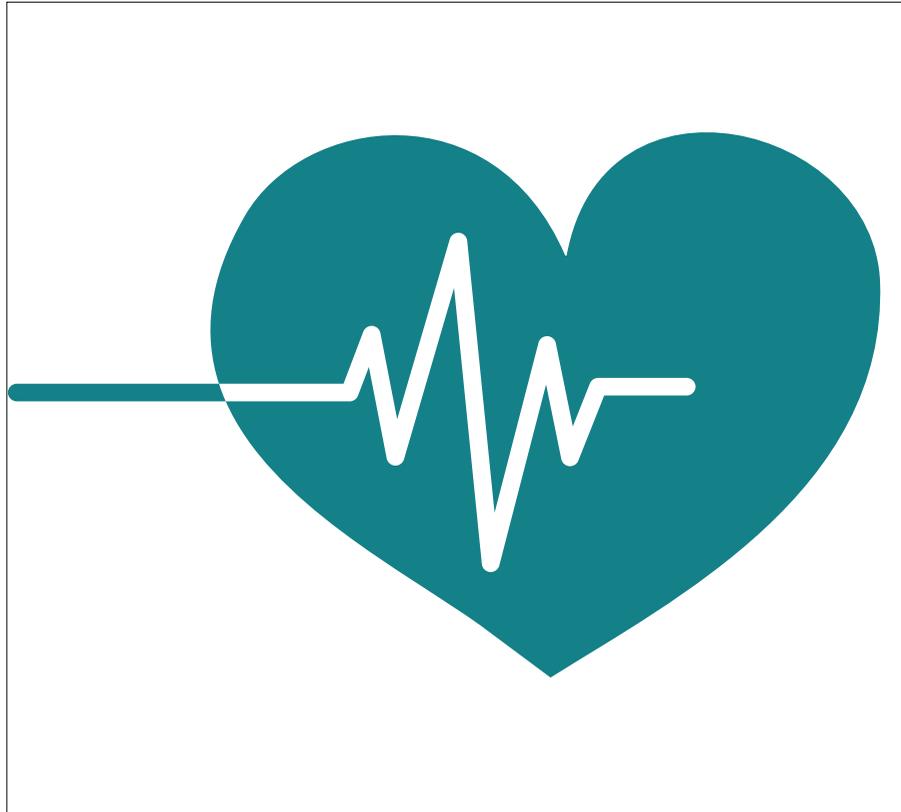
HTRF solutions for cardiology

Revvity's kits and services offer simple, rapid, cost-effective solutions for cardiology research.

Referenced in a growing number of publications aiming at understanding the pathophysiology of atherosclerosis, heart fibrosis, and cardio protection.

Comprehensive portfolio of Total and Phospho-protein, cytokine, kinase and biomarker assays, covering all major pro-pathways, including Nitric Oxide signaling.

Large selection of GPCR tools and solutions, including our acclaimed **cAMP** and **IP-One** assays.



GPCRs	
<i>cAMP – Gs Dynamic</i>	<i>Tag-lite beta1</i>
<i>cAMP – Gs HiRange</i>	<i>Tag-lite beta2</i>
<i>IP-One – Gq</i>	<i>Tag-lite AT2</i>
<i>Forskolin</i>	<i>Tag-lite A2A</i>
<i>IBMX</i>	<i>Tag-lite A2B</i>
	<i>Tag-lite A1</i>
	<i>Tag-lite A3</i>

TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)

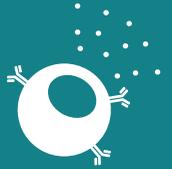
DISEASES	Hypertension Hypercholesterol Atherosclerosis Acute and Chronic Ischemia	Atherosclerosis	Stroke Hypertrophy	Heart fibrosis
	NITRIC OXIDE	APOLIPOPROTEINSE	CARDIO PROTECTION	FIBROSIS
Cytokines 	<i>IFN gamma</i> <i>TNF alpha</i> <i>IL2</i> <i>IL4</i> <i>IL5</i> <i>IL6</i> <i>IL10</i> <i>IL12</i> <i>IL17</i>			<i>TGF beta</i> <i>IL6</i> <i>IL1 beta</i> <i>TNF alpha</i>
Phospho-proteins 	<i>Alpha Tubulin</i> <i>Phospho-VASP (Ser157)</i> <i>Phospho-VASP (Ser239)</i> <i>Total VASP</i> <i>Total & Phospho-Phospholamban (Thr17)</i>		<i>Alpha Tubulin</i> <i>Phospho-IRS1 (Ser307 mouse/Ser312 human)</i> <i>Total & Phospho-AKT1 (Ser473)</i> <i>Total & Phospho-AKT2 (Ser473)</i> <i>Total & Phospho-AKT3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Thr308)</i> <i>Total AKT1/2/3</i> <i>Phospho-mTOR (Ser2448)</i> <i>Total & Phospho-GSK3 alpha (Ser21)</i> <i>Phospho-BAD (Ser112)</i> <i>Total & Phospho-ERK (Thr202/Tyr204)</i>	<i>Alpha Tubulin</i> <i>Total & Phospho-SMAD1 (Ser463/465)</i> <i>Phospho-SMAD2 (Ser465/427)</i> <i>Total & Phospho-SMAD3 (Ser423/425)</i>
Biomarkers 	<i>cGMP</i>	<i>PCSK9</i> <i>APO-E/A1/B/C3</i>		<i>HA</i> <i>P3NP</i> <i>Pro-collagen</i> <i>MMP2/9 & TIMP-1</i>

TABLE OF CONTENTS

HTRF solutions for autoimmunity

HTRF solutions for cardiology

HTRF solutions for NAFLD and fibrosis

HTRF solutions for neuroscience

HTRF solutions for CAR-T research

HTRF solutions for TLR research

HTRF solutions for virology

NAFLD AND FIBROSIS

HTRF solutions for NAFLD and fibrosis

Revvity's kits and services offer simple, rapid, cost-effective solutions for NAFLD and fibrosis.

Successfully applied to all the steps preceding and following the state of non-alcoholic steatohepatitis.

Comprehensive portfolio of:

- Cytokine assays
- Chemokine assays
- Kinase assays

covering all major pro-inflammatory pathways. Of particular interest are our star products in fibrosis, including **Alpha-SMA**, **Pro-collagen**, and **Hyaluronic acid**.

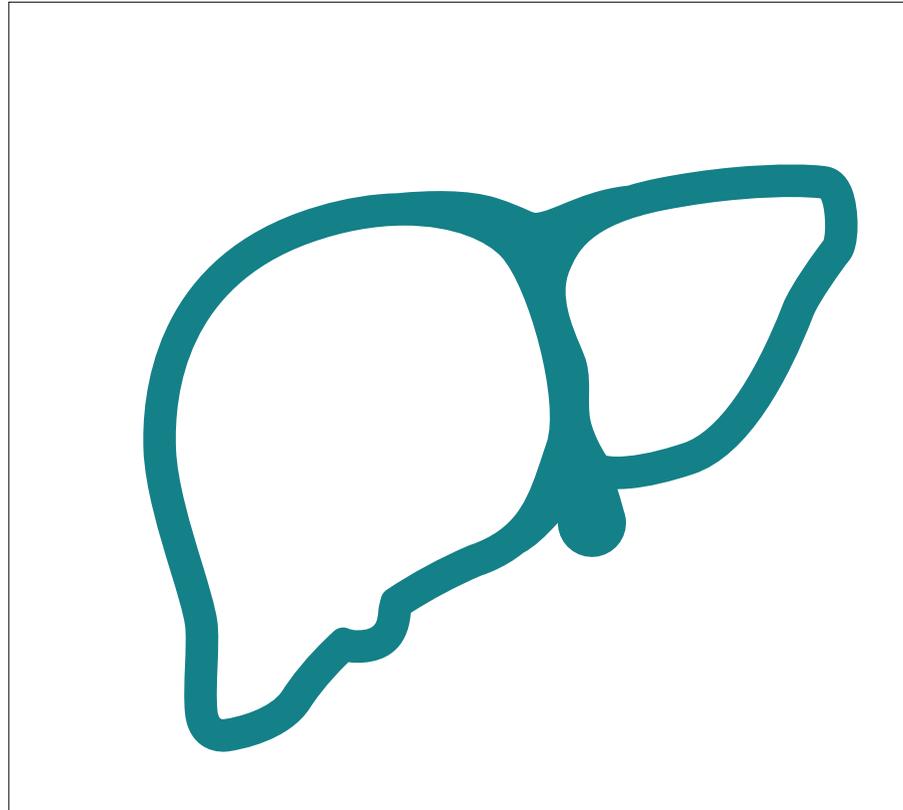


TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)

	NUCLEAR RECEPTORS	STEATOSIS	IR and DNL	ER STRESS	FIBROSIS
 Cytokines			IL1 beta IL6 TNF alpha IFN gamma		TGF beta CCL2 IL2 IL18
 Phospho-proteins		Alpha Tubulin Total & Phospho-FoxO1 (Ser256) Total & Phospho-ACC (Ser79) Total & Phospho-AMPK (Trh172) Total & Phospho-CREB (Ser133)	Alpha Tubulin Total & Phospho-FoxO1 (Ser256) Total & Phospho-ACC (Ser79) Total & Phospho-AMPK (Trh172) Total & Phospho-CREB (Ser133) Total & Phospho-IRS1 (Ser307) Total & Phospho-SMAD3 (Ser423/425) Total & Phospho-IR beta (Tyr1150/1151) Phospho-4EBP1 (Thr37/48) Total & Phospho-GSK3 beta (Ser9)	Alpha Tubulin Total & Phospho-EIF2 alpha (Ser52) Total & Phospho-FoxO1 (Ser256)	Alpha Tubulin Total & Phospho-SMAD3 (Ser423/425) Total & Phospho-SMAD2 (Ser465/427) Phospho-PDGFRβ beta (Tyr751) Phospho-P38 (Thr180/Tyr182) Total & Phospho-NF-kB (Ser536) Total & Phospho YAP (Ser127) Total & Phospho-FAK (Tyr397)
 Biomarkers		ApoC3,E,A1,B	Insulin High range Insulin Ultra sensitive Mouse Insulin serum		Hyaluronic acid Pro-collagen Alpha-SMA MMP2,9/TIMP-1 P3NP
 PPis	Streptavidin-XL665 MAb Anti FLAG M2-d2 MAb Anti FLAG M2-Eu cryptate MAb Anti FLAG M2-Tb cryptate MAb Anti FLAG M2-XL665 MAb Anti-6HIS-Eu cryptate Gold MAb Anti-6HIS Tb cryptate Gold MAb Anti 6HIS-d2 MAb Anti 6HIS-XL665 MAb Anti GST-d2 MAb Anti GST-Eu cryptate MAb Anti GST-Tb cryptate MAb Anti GST-XL665				

TABLE OF CONTENTS

HTRF solutions for autoimmunity

HTRF solutions for cardiology

HTRF solutions for NAFLD and fibrosis

HTRF solutions for neuroscience

HTRF solutions for CAR-T research

HTRF solutions for TLR research

HTRF solutions for virology

NEUROSCIENCE

HTRF solutions for neuroscience

Revvity's kits, including total and Phospho-proteins, Cytokines, and Biomarkers, offer simple, rapid, cost-effective solutions to researchers working in neurodegenerative diseases.

A comprehensive product pipeline and a growing number of publications aiming at understanding the pathophysiology of Alzheimer's, Parkinson's, ALS/FTD diseases onset.

A full range of HTRF assays to study chronic neuroinflammation triggered by glial cells and delineate their critical role in neurodegeneration pathways.



TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)

	ALZHEIMER'S	PARKINSON'S	ALS/FTD	MICROGLIA	ASTROCYTES
 Cytokines	IL1β IL4 IL6 IL10 TNF alpha	IL1β IL6 TNF alpha	IL1β IL6 TNF alpha IL8	IL1β IL4 IL6 IL10 TGF betaβ IL2 IL18	IL1β IL10 TNF alpha IL17 CXCL1 CCL2 TGFβ beta
 Phospho-proteins	Alpha Tubulin Total & Phospho-GSK3 βbeta (Ser9) Total & -Phospho Tau (Ser202) Phospho & Total-Tau (Ser404) Phospho-AKT1/2/3 (Ser473) Phospho-AKT1/2/3 (Thr308) Total AKT1/2/3 Total & Phospho-ERK (Thr202/Tyr204)	Alpha Tubulin Total & Phospho-aSyn (Ser129) Total & Phospho-LRRK2 (Ser935) Phospho-mTOR (Ser2448) Phospho-ULK1 (Ser757/Ser758) Total & Phospho-AMPK (Trh172) Total & Phospho-TBK1 (Ser172) Phospho-BAD (Ser112) Phospho-P38 (Thr180/Tyr182) Phospho-JNK (Thr183/Tyr185) Total & Phospho-MKK4 (Ser257) Phospho-AKT1/2/3 (Ser473) Phospho-AKT1/2/3 (Thr308) Total AKT1/2/3 Total & Phospho-ERK (Thr202/Tyr204) Total & Phospho-P62SQSTM1 (Ser403)	Alpha Tubulin Phospho-mTOR (Ser2448) Phospho-ULK1 (Ser757/Ser758) Total & Phospho-AMPK (Trh172) Total & Phospho-TBK1 (Ser172) Phospho-BAD (Ser112) Total & Phos- pho-P62SQSTM1 (Ser403)	Alpha Tubulin Total & Phospho-SLP76 (Ser376) Total & Phospho-BTK (Tyr223) Phospho-AKT1/2/3 (Ser473) Phospho-AKT1/2/3 (Thr308) Total AKT1/2/3 Total & Phospho-ERK (Thr202/Tyr204) Total & Phospho-MKK4 (Ser257) Total & Phospho-NF-kB (Ser536)	Alpha Tubulin Total & Phospho-IKKbeta (Ser177/181) Total & Phospho-NF-kB (Ser536) Total & Phospho-MEK1 (Ser298) Total & Phospho-MEK1/2 (Ser218/222) Total & Phospho-ERK (Thr202/Tyr204) Total & Phospho-STAT3 (Tyr705)
 Biomarkers	Tau aggregation MAP-TAU Amyloid beta1-40	a-Synuclein aggregation BDNF Cleaved PARP	Cleaved PARP Progranulin	TREM2 HMGB1	MMP9 MANF BDNF
 PPIs/ Tag-lite/ GPCRs	Biotin-APP-Eu cryptate APP	Tag-Lite D1 Receptor Tag-Lite D2 Receptor		IP-one	IP-one

TABLE OF CONTENTS

HTRF solutions for autoimmunity

HTRF solutions for cardiology

HTRF solutions for NAFLD and fibrosis

HTRF solutions for neuroscience

HTRF solutions for CAR-T research

HTRF solutions for TLR research

HTRF solutions for virology

CAR-T RESEARCH

HTRF solutions for CAR-T research

Revvity's Total and Phospho-protein, Cytokine, and biomarker kits and our services offer simple, rapid, cost-effective solutions to researchers working in immuno-oncology.

Early drug discovery stages on CAR-T engineering and functional testings can be addressed with our full range of HTRF assays.

Early DD

- Address tumor targeting
- Intrinsic T cell potency: Functionality and persistence
- Extrinsic T cell potency: Action on the tumor microenvironment

INTRINSIC T CELL POTENCY

	Assess functionality	Assess persistence
 Cytokines	<i>IFN gamma</i> <i>TNF alpha</i> <i>IL1 beta</i> <i>IL2</i> <i>IL6</i> <i>IL12</i> <i>IL18</i> <i>GM-CSF</i>	
 Phospho-proteins	<i>Alpha Tubulin</i> <i>Total & Phospho-ZAP70 (Tyr319)</i> <i>Total & Phospho-SLP76 (Ser376)</i> <i>Total & Phospho-SHP1 (Tyr564)</i> <i>Total & Phospho-SHP2 (Tyr542)</i> <i>Total & Phospho-STAT1 (Tyr701)</i> <i>Phospho-STAT5 (Tyr694)</i> <i>Phospho-STAT6 (Tyr641)</i> <i>Total & Phospho-ERK (Thr202/Tyr204)</i> <i>Total & Phospho-AKT1 (Ser473)</i> <i>Total & Phospho-AKT2 (Ser473)</i> <i>Total & Phospho-AKT3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Ser473)</i> <i>Phospho-AKT1/2/3 (Thr308)</i> <i>Total AKT1/2/3</i> <i>Phospho mTor (Ser2448)</i> <i>Phospho JNK (Thr183/Tyr185)</i> <i>Phospho-cJun (Ser63)</i> <i>Total & Phospho-NF-kB (Ser536)</i> <i>Phospho/Total-CREB (Ser133)</i>	
 Biomarkers	<i>Ki67</i> <i>GRZB</i> <i>Perforin</i>	<i>Ki67</i> <i>BAD</i> <i>Cleaved PARP</i>

TABLE OF CONTENTS

HTRF solutions for autoimmunity

HTRF solutions for cardiology

HTRF solutions for NAFLD and fibrosis

HTRF solutions for neuroscience

HTRF solutions for CAR-T research

HTRF solutions for TLR research

HTRF solutions for virology

TLR RESEARCH

HTRF solutions for TLR research

Revvity's Cytokine, Total and Phospho-protein kits as well as services offer simple, rapid, cost-effective solutions to researchers in inflammatory related diseases.

Drugs targeting innate immunity especially Toll-like receptor signaling can be addressed with our full range of HTRF assays.

	TLR signaling
 <p>Cytokines</p>	<p><i>TNF alpha</i> <i>IL6</i> <i>IL1 beta</i> <i>IFN alpha</i> <i>IFN beta</i></p>
 <p>Phospho-proteins</p>	<p><i>Alpha Tubulin</i> <i>Total IRAK3</i> <i>Total & Phospho-TBK1 (Ser172)</i> <i>Total & Phospho-IKKbeta (Ser177/181)</i> <i>Total & Phospho-NF-kB (Ser536)</i> <i>Phospho-IRF3 (Ser386)</i> <i>Phospho-JNK (Thr183/Tyr185)</i> <i>Phospho-cJun (Ser63)</i> <i>Phospho-P38 (Thr180/Tyr182)</i></p>

TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)

VIROLOGY

HTRF solutions for virology

Revvity's PPI reagents, kits, and services offer simple, rapid, cost-effective solutions to researchers working in virology.

They have been successfully applied to the discovery of anti-viral agents targeting the binding, integration, replication, and assembly steps of the viral cycle.

Renowned companies have published on their use of Revvity products to develop creative approaches to antiviral drug development. Among them are Medimmune, the Scripps Institute, and the NIH.

HTRF reagents have been used to look for inhibitors of HRV attachment and replication, and inhibitors of HIV capsid formation and integration.

	Other products of interest
Custom services	<p><i>Custom labelling</i></p> <p><i>Custom assay development</i></p>
Specialty assays	<p><i>Viral double-stranded RNA detection kit</i></p>

PPI reagents
<p><i>MAb Anti HA-d2</i></p> <p><i>MAb Anti HA-Eu cryptate</i></p> <p><i>MAb Anti HA-Tb cryptate</i></p> <p><i>MAb Anti HA-XL665</i></p>
<p><i>MAb Anti DNP-Eu cryptate</i></p> <p><i>MAb Anti DNP-XL665</i></p>
<p><i>MAb Anti FLAG M2-d2</i></p> <p><i>MAb Anti FLAG M2-Eu cryptate</i></p> <p><i>MAb Anti FLAG M2-Tb cryptate</i></p> <p><i>MAb Anti FLAG M2-XL665</i></p>
<p><i>MAb Anti GST-d2</i></p> <p><i>MAb Anti GST-Eu cryptate</i></p> <p><i>MAb Anti GST-Tb cryptate</i></p> <p><i>MAb Anti GST-XL665</i></p>
<p><i>MAb Anti-6HIS-Eu cryptate Gold</i></p> <p><i>MAb Anti-6HIS Tb cryptate Gold</i></p> <p><i>MAb Anti 6HIS-d2</i></p> <p><i>MAb Anti 6HIS-XL665</i></p>
<p><i>MAb Anti cmyc-d2</i></p> <p><i>MAb Anti cmyc-Eu cryptate</i></p> <p><i>MAb Anti cmyc-Tb cryptate</i></p> <p><i>MAb Anti cmyc-XL665</i></p>

TABLE OF CONTENTS

[HTRF solutions for autoimmunity](#)

[HTRF solutions for cardiology](#)

[HTRF solutions for NAFLD and fibrosis](#)

[HTRF solutions for neuroscience](#)

[HTRF solutions for CAR-T research](#)

[HTRF solutions for TLR research](#)

[HTRF solutions for virology](#)



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.

317157