

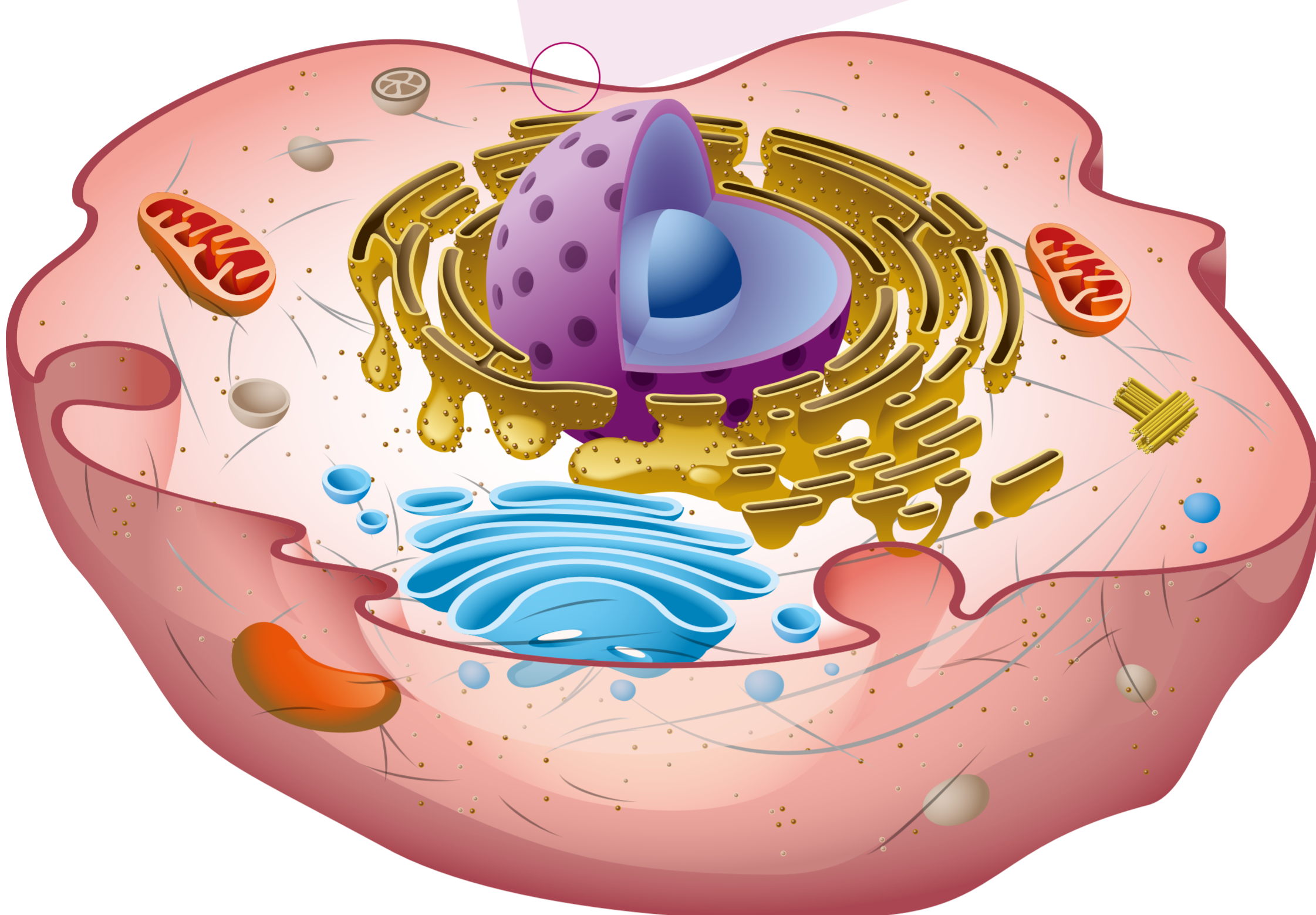
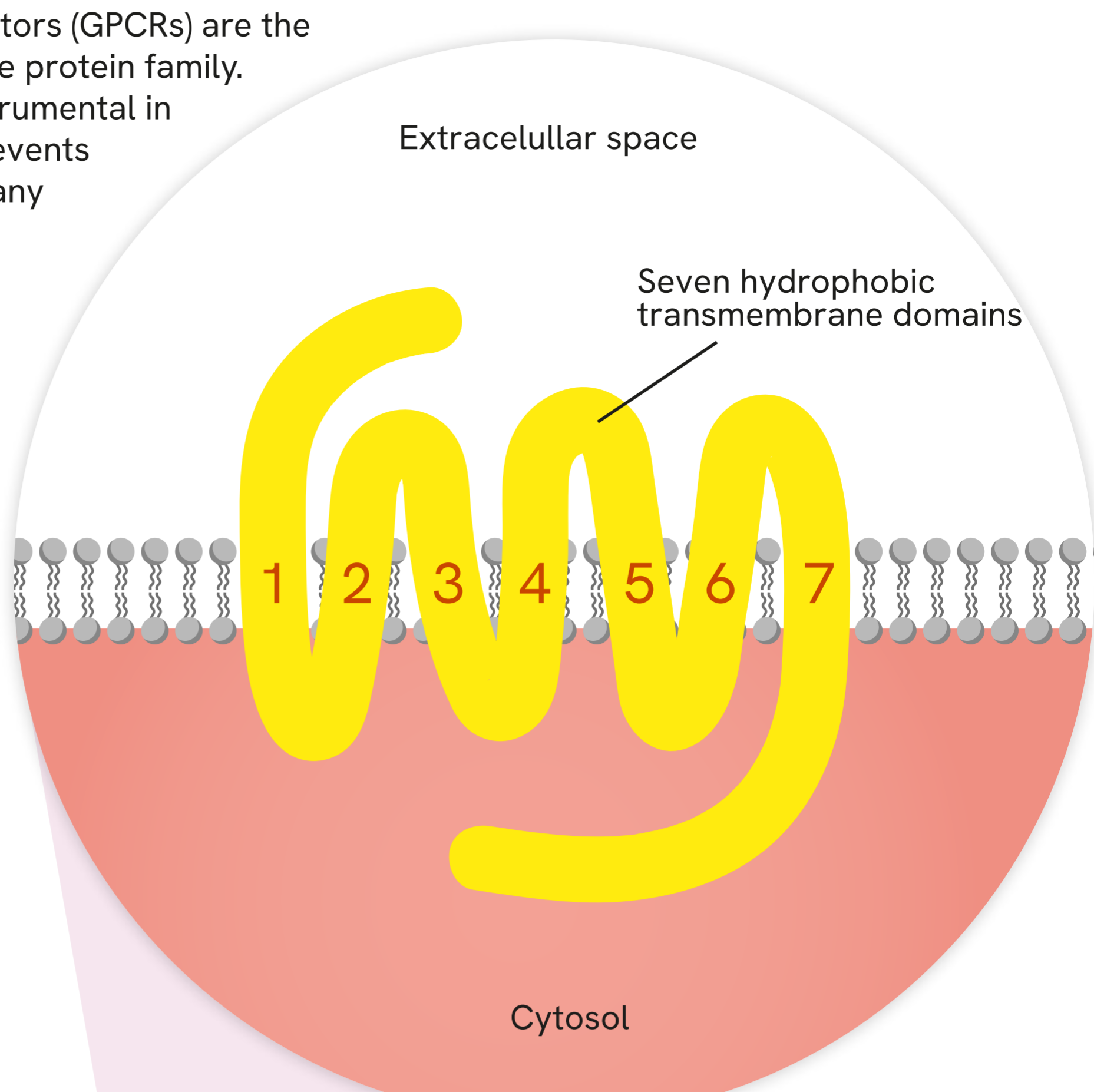
GPCRs and their prevalence in drug therapies.

G protein-coupled receptors (GPCRs) are the largest human membrane protein family. These receptors are instrumental in numerous cell signaling events and play a key role in many physiological processes and pathological conditions.^{1,2}

GPCRs can associate with a wide variety of ligands including...

- Hormones
- Neurotransmitters
- Lipids
- Nucleotides
- Ions

GPCR signaling pathways can be activated by canonical ligands or biased ligands



GPCR signaling via G-proteins

1. In the presence of a canonical ligand the inactive GPCR is "stimulated" altering its conformation and its association with the alpha subunit ($G\alpha$) of the G-protein.

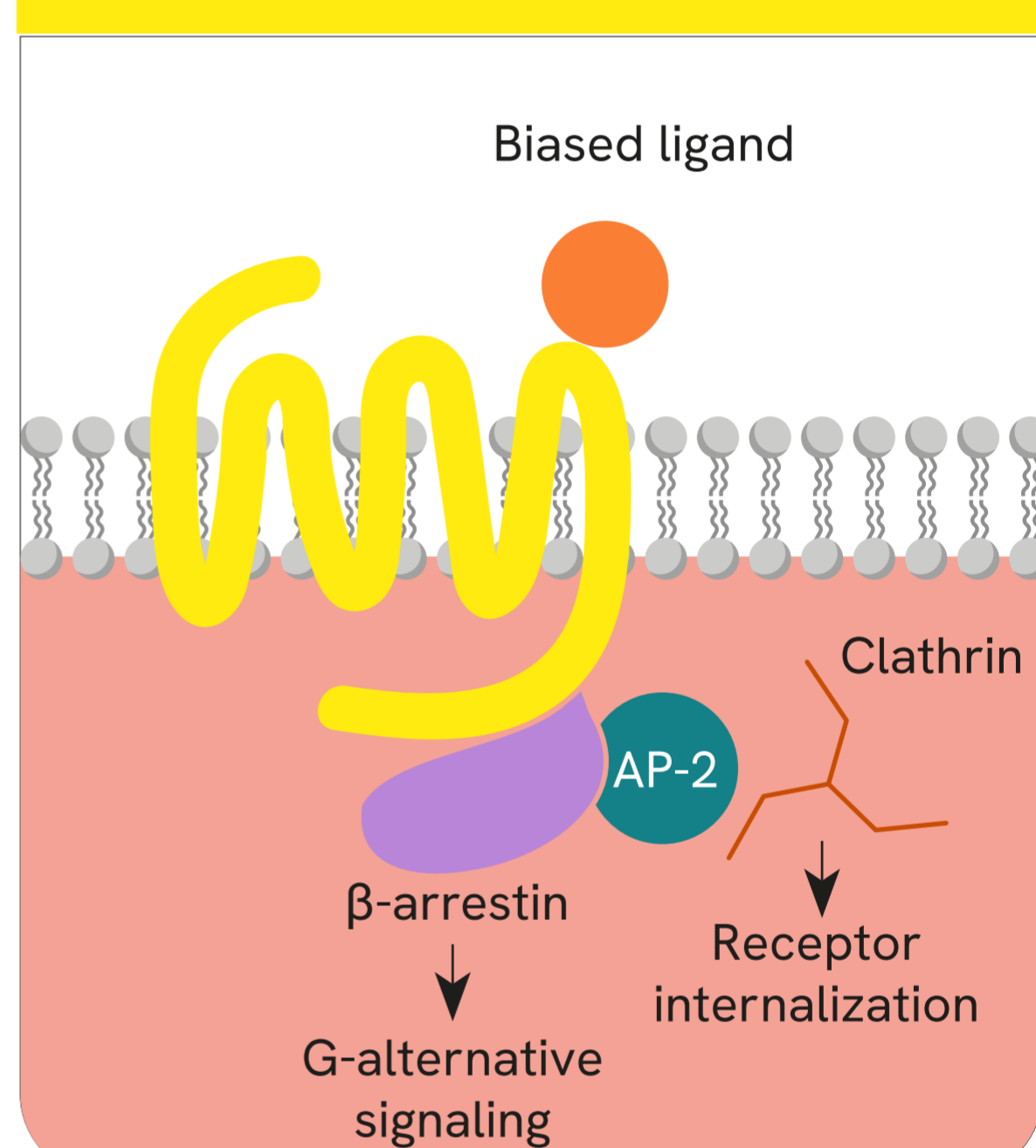
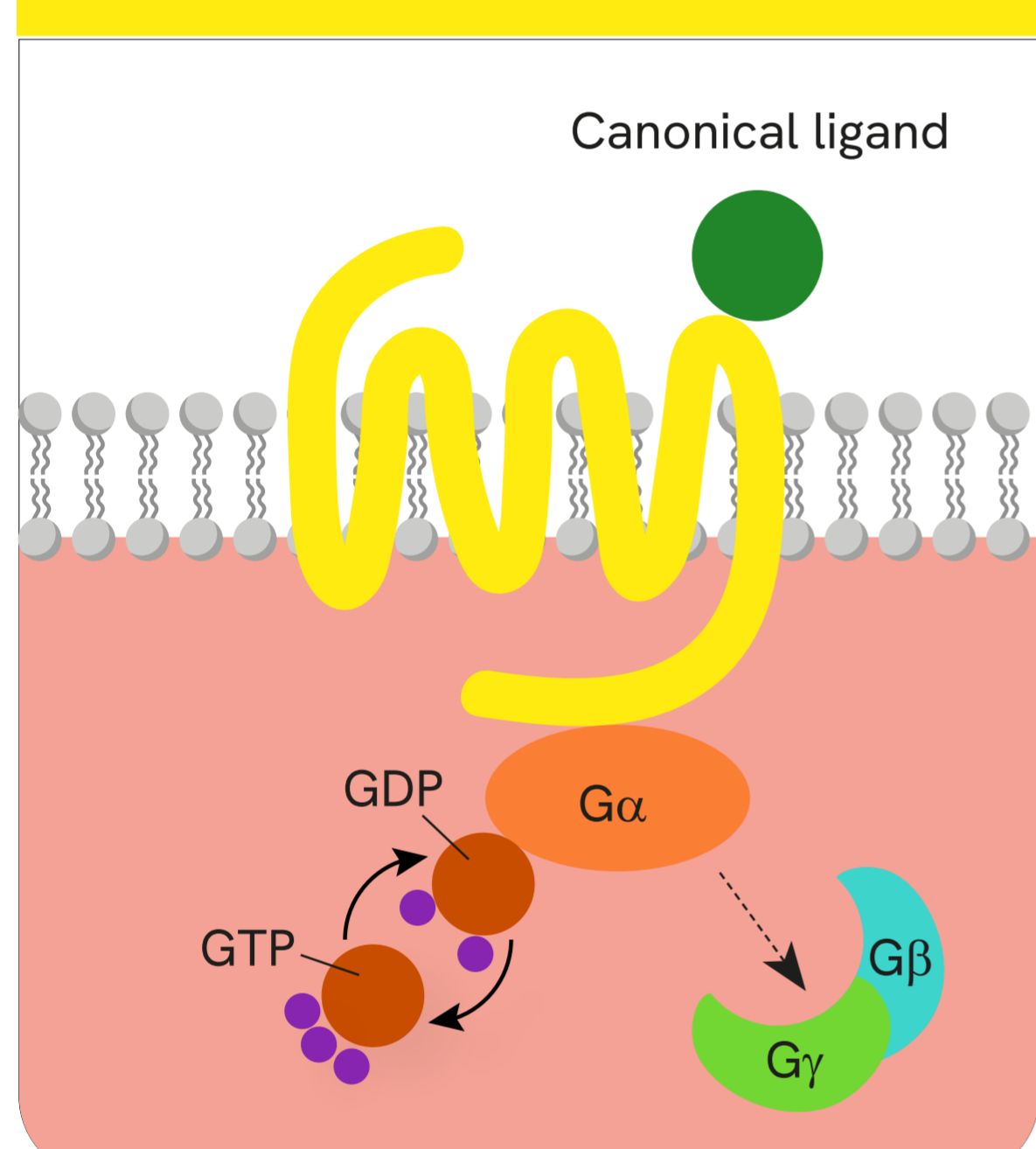
2. GDP is exchanged for GTP and $G\alpha$ dissociates from the beta ($G\beta$) and gamma ($G\gamma$) subunits.

3. The dissociated $G\alpha$ and $G\beta$ - $G\gamma$ dimer can now interact with other proteins, stimulating cell signaling pathways.

Arrestin-mediated signaling

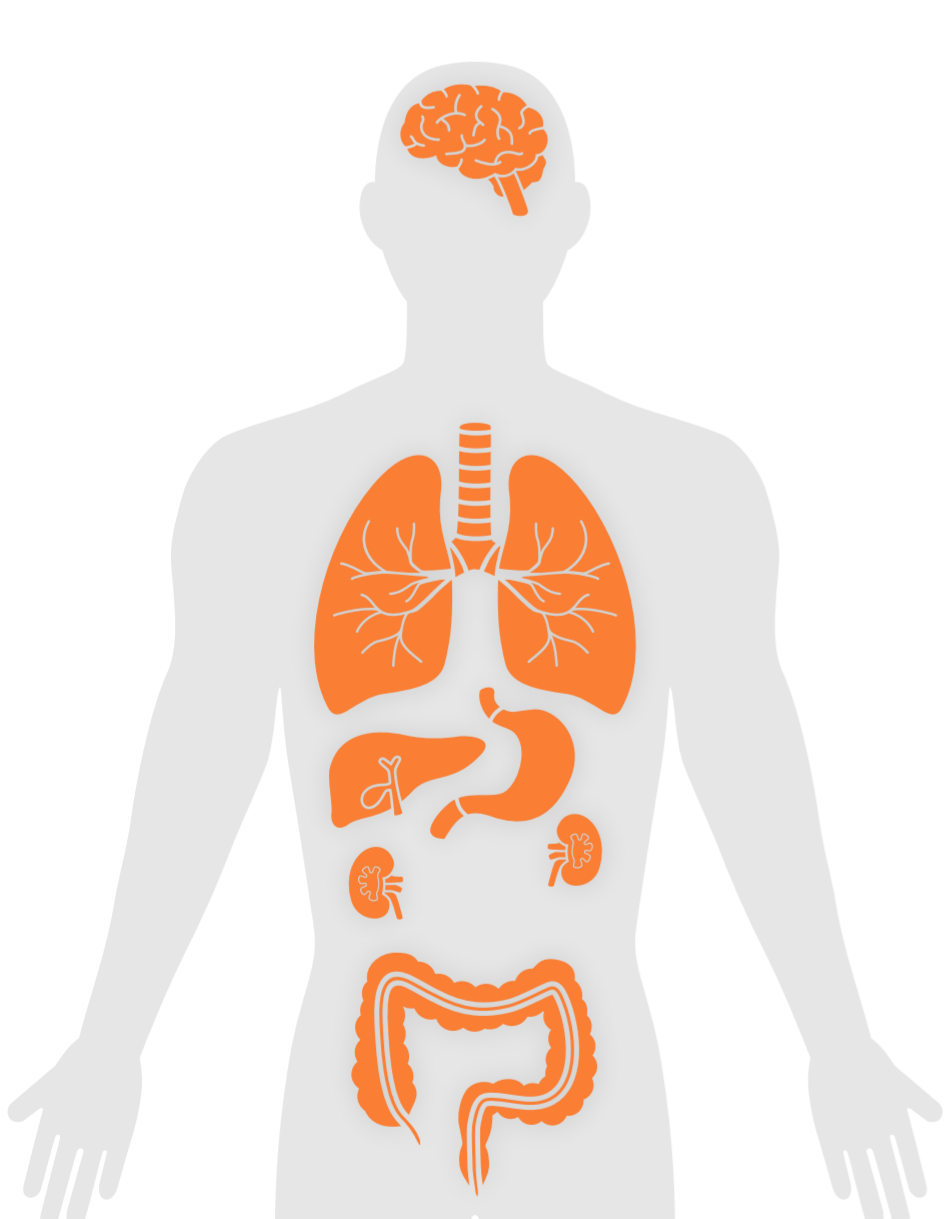
Arrestins can influence GPCR signaling in several ways.

In addition to competitively blocking the binding of G-proteins, arrestins also regulate GPCRs by facilitating their endocytosis via internalization into clathrin-coated pits.³



Most of the body's GPCRs are controlled by just two arrestin proteins, β -arrestin 1 and β -arrestin 2.

The flexible structure and versatile functions of arrestins make them highly pliable drug targets which, in time, will help to fine tune the disease-specificity of GPCR drugs, making them safer and more effective.



GPCR-targeted drugs already exist for numerous indications.^{4,5,6,7}

Cardiovascular diseases	Atropine (Bradycardia treatment) <i>Competitive antagonist of acetylcholine muscarinic receptors</i>
Neuroscience	Carvedilol (Congestive heart failure/hypertension treatment) <i>Inverse agonist of the β-adrenergic receptor</i> Bupropion (Nicotine withdrawal treatment) <i>Partial agonist of acetylcholine nicotinic receptors</i> Morphine (Pain treatment) <i>Full agonist of μ-opioid receptors</i>
Metabolic diseases	Exenatide (Type 2 diabetes treatment) <i>Agonist of the glucagon-like peptide-1 receptor</i>
Oncology	Degarelix (Prostate cancer treatment) <i>Antagonist of gonadotropin-releasing hormone</i> Vismodegib (Basal cell carcinoma treatment) <i>Competitive antagonist of Smoothed (Smo)</i>
Immunology	CF 101 (Rheumatoid arthritis treatment) <i>Antagonist of the adenosine A3 receptor</i>

1. Hauser AS, Attwood MM, Rask-Andersen M, Schiöth HB, & Gloriam DE. Trends in GPCR drug discovery: new agents, targets and indications. *Nature Reviews Drug Discovery*. 2017;16(12):829-842. doi: 10.1038/nrd.2017.178 2.
2. Essentials of cell biology: g-protein-coupled receptors play many different roles in eukaryotic cell signaling. *Nature Education*. <https://www.nature.com/scitable/e-books/essentials-of-cell-biology-14749010/122997540>. Published 2014. Accessed May 21, 2021.
3. Mj L, C H. Arrestin interactions with G protein-coupled receptors. *Handb Exp Pharmacol*. 2014;219:15-56. doi: 10.1007/978-3-642-41199-1_

4. Pharmacology: Insight Into Ligand Classes. Revvity. Accessed May 21, 2021.
5. Exenatide. BNF: British National Formulary - NICE. <https://bnf.nice.org.uk/drug/exenatide.html#drugAction>. Accessed May 23, 2021.
6. Aditya S, Rattan A. Vismodegib: A smoothed inhibitor for the treatment of advanced basal cell carcinoma. *Indian Dermatol Online J*. 2013;4(4):365-368. doi: 10.4103/2229-5178.120685
7. Plicidenoson. DrugBank Online. <https://go.drugbank.com/drugs/DB055111>. Published 2007. Updated February 21, 2021. Accessed May 23, 2021.