

GPCR Efficacy and Biased Agonism

Patrick M. Sexton^{1,2}, Arthur Christopoulos¹, Denise Wootten^{1,2}

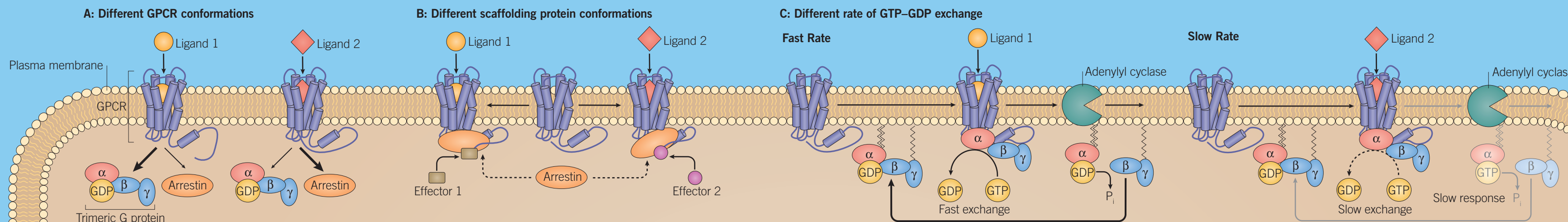
¹Drug Discovery Biology, Monash Institute of Pharmaceutical Sciences, Monash University, Parkville 3052, Victoria, Australia

²School of Pharmacy, Fudan University, Shanghai 2012013, China.

TOCRIS
a biotechne brand

Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals

www.tocris.com

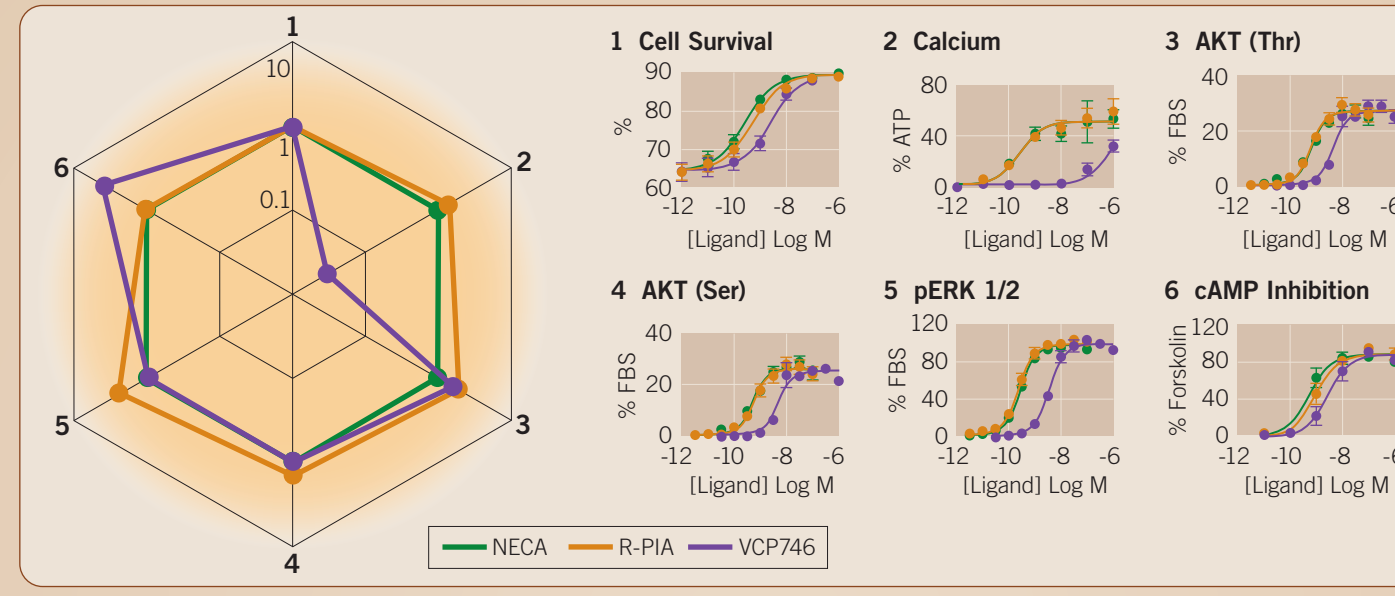


Multiple Mechanisms Elicit Biased Agonism

(A) Different ligands induce different conformations within G protein-coupled receptors (GPCRs), promoting binding conditions for particular transducer proteins such as G proteins and arrestins. (B) Specific ligand-induced receptor conformations induce different conformational changes within scaffolding proteins such as arrestins, causing activation of distinct downstream signaling pathways. (C) Different ligands can control the rate of GTP-GDP exchange through induction of distinct conformational rearrangements within G proteins. Ligands that promote fast GTP association and G protein turnover allow more signaling events to occur than ligands that induce a slow rate of exchange.

G Protein-Coupled Receptors (GPCRs)

GPCRs are seven transmembrane domain receptors, which can move between inactive-like (R , R' and R'') and active-like (R^* and R^{**}) states. This may occur in the absence of ligand or transducer (the Apo state), however the energy barrier to achieving and maintaining active-like conformations means that it's unlikely to occur or has a lower probability in the absence of an agonist and G protein (or other transducer). The presence of both an agonist and transducer (e.g. G protein) stabilizes the active GPCR conformation and gives the highest probability that a signaling event will occur. The rates of transducer engagement and turnover are altered in an agonist dependent manner that determines observed efficacy¹. Moreover, GPCRs can interact with multiple distinct transducers or regulatory proteins and these can be preferentially engaged in an agonist-specific manner giving rise to biased agonism¹. Nanobodies that bind to the intracellular face of GPCRs can allosterically drive either inactive or active-like conformations and have been used in structural studies to capture these states², however increasing numbers of structures with G proteins are now being solved³.

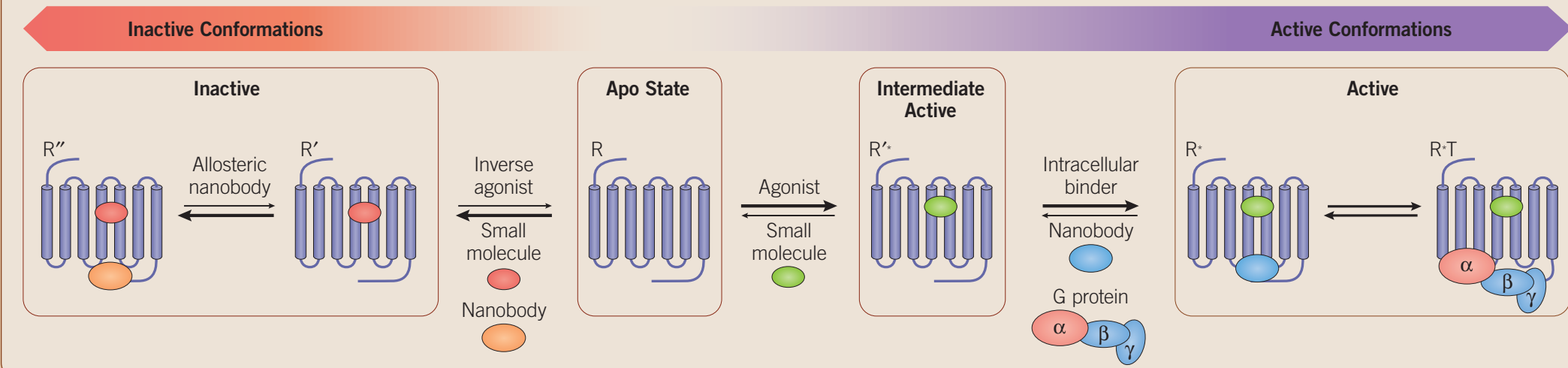


Identifying Biased Agonism at the Adenosine A₁ GPCR (A₁R) Using the Operational Model

The six concentration-response curves (pictured) show activation of different signaling events for three A₁R agonists, NECA (green line), R-PIA (orange line) and VCP746 (purple line)⁴. Operational modeling was used to derive the transduction ratio ($\log\tau/K_A$; linking agonist occupied receptors to pharmacological effects) and this was plotted in a "Web of Bias"⁵.

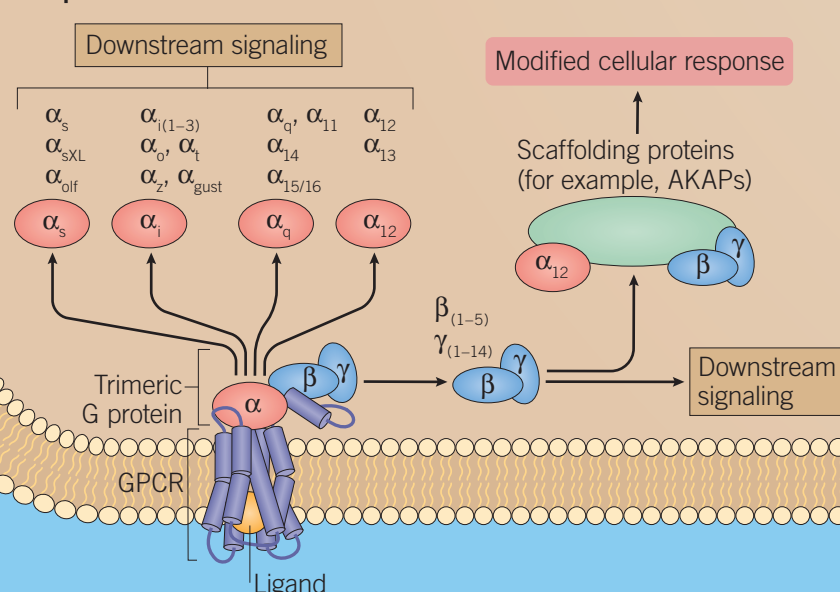
The "Web of Bias" chart shows each drug profile relative to the reference agonist NECA and reference pathway 1 (as $\Delta\Delta\log\tau/K_A$). Using this visual representation, the biased agonism of VCP746 for pathway 2 (reduced response relative to NECA and pathway 1), is easily observed. In contrast, R-PIA displays a similar signaling profile to the reference agonist, NECA.

Of note, biased agonism is always relative, and must be expressed by comparison with a reference ligand; observations of bias can differ in a system (cell background)-dependent, and temporal manner¹.

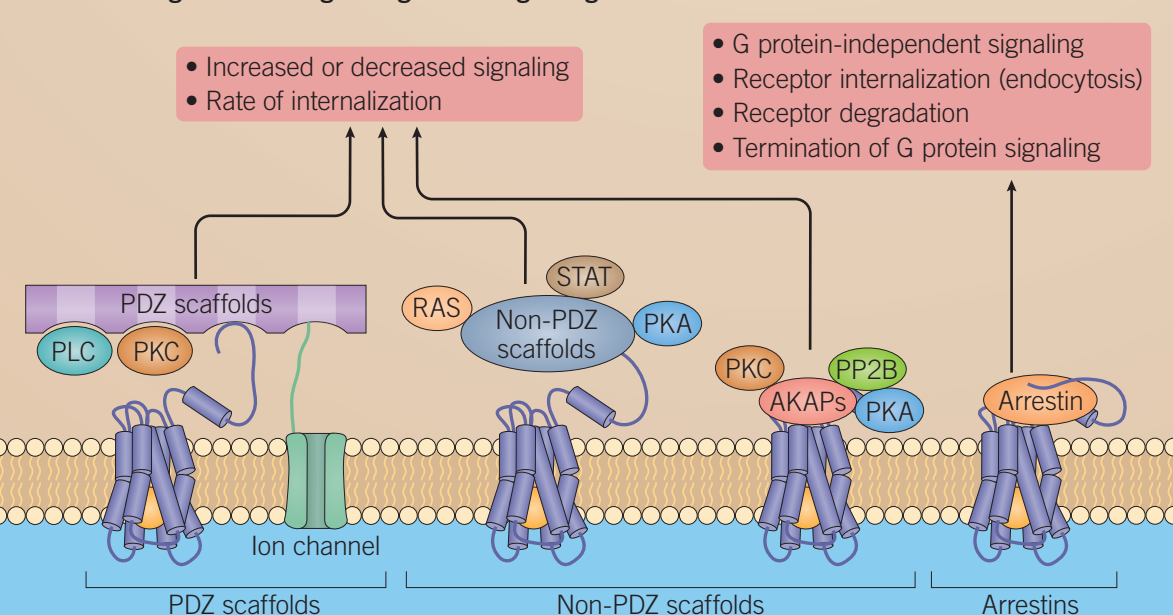


GPCRs are pleiotropically coupled and so can partner with multiple different G proteins (A) or other protein transducers (B) to regulate intracellular signaling.

A: G protein Activation

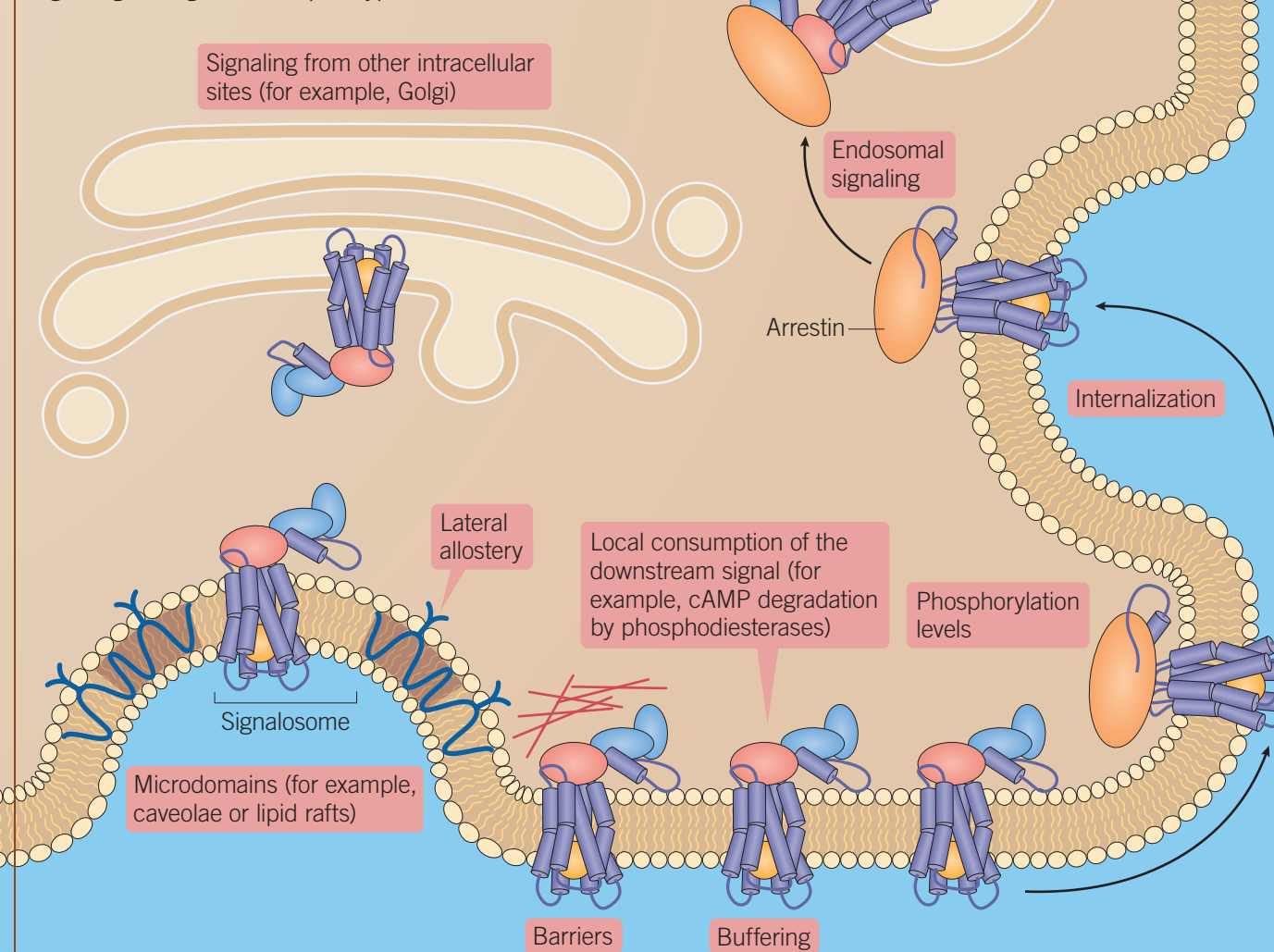


B: Scaffolding Proteins Regulating GPCR Signaling



Spatial and Kinetic Factors

Biased agonism is regulated by the temporal stability of biomolecular interactions between ligand, receptor and transducers or regulatory proteins. Through biased agonism, different ligands can alter the location, type and strength of signaling through one receptor type.



Products available from Tocris

- Key**
- Agonist
 - Antagonist
 - Inverse Agonist
 - Partial Agonist
 - Positive Allosteric Modulator
 - Negative Allosteric Modulator

- 5-HT Receptors**
- WAY 181187
 - SB 269970
 - SB 243213
 - Tansospirone
 - PU 02

- Adenosine A₁ Receptors**
- (+)-5'-Chloro-5'-deoxy-ENBA
 - DPCPX
 - PQ 69
 - PD 81723

- Adenosine A_{2A} Receptors**
- CGS 21680
 - Istradefylline
 - LUF 5834

- Adenosine A_{2B} Receptors**
- BAY 60-6583
 - PSB 603
 - MRS 1706
 - LUF 5834

- Adenosine A₃ Receptors**
- 2-Cl-IB-MECA
 - MRS 1220
 - PSB10

- Adrenergic Receptors**
- A 61603
 - SR 59230A
 - ICI 118,551
 - Xamoterol

- Cannabinoid Receptors**
- HU 308
 - AM 251
 - AM 630
 - Cannabigerol
 - Org 27569
 - PSNCBAM-1

- Dopamine Receptors**
- SKF 81297
 - L-741,626
 - Fenoldopam
 - PAOPA

- GABA Receptors**
- (R)-Baclofen
 - CGP 55845
 - MRK 016
 - TP 003
 - CGP 7930

- Glutamate (Metabotropic) Receptors**
- (S)-3,5-DHPG
 - LY 341495 disodium salt
 - VU 0360172
 - ADX 10059

- Muscarinic Receptors**
- Xanomeline
 - (S)-(+)-Dimethindene
 - VU 0238429

- Opioid Receptors**
- SNC 80
 - Naltrindole
 - Meptazinol
 - BMS 986187

- Purinergic Receptors**
- ATPyS tetralithium salt
 - MRS 2500
 - GW 791343

- Trace Amine 1 Receptor**
- Ractopamine
 - EPPTB

- Heterotrimeric G-protein GTPase**
- SCH 202676
 - NF 023
 - G-Protein antagonist peptide
 - CMPD101

- DREADDs**
- DREADD agonist 21
 - Perlapine
 - Salvinorin B
 - Clozapine N-oxide dihydrochloride

- References**
- Wootten et al. (2018) *Nat Rev Mol Cell Biol* 19:638-653
 - Thal et al. (2018) *Nature* 559:45-53
 - Glukhova et al. (2018) *ACS Pharmacol Transl Sci* 1:73-83
 - Baltos et al. (2016) *Biochem Pharmacol* 99:101-112
 - Kenakin et al. (2012) *ACS Chem Neurosci* 3:193-203

Therapeutic Application of Biased Agonism

Studying the pharmacology of GPCR biased agonists can help improve the efficacy and safety profiles of drugs. Identification of compounds with a therapeutically beneficial activation profile that have reduced undesired on-target signaling can improve patient outcomes and translation of drugs to the clinic. However, the selection of the most appropriate biased agonists will require improved understanding of

the patterns of tissue-specific and integrated signaling that mediate different outcomes both physiologically and in a disease specific context. Collectively, such advances in understanding of disease, combined with advances in drug design, GPCR structure determination and pharmacological assessment of compounds should provide increasing numbers of new biased agonist drugs to the clinic.

For copies of this poster, please visit tocris.com
© 2019 Tocris Cookson, Ltd. Tocris is a Bio-Techne brand