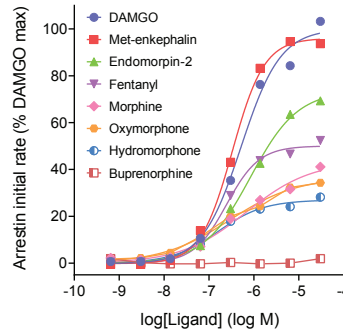




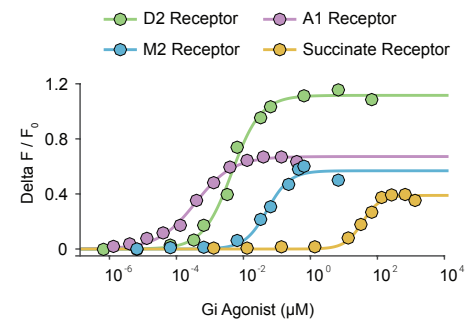
Montana Molecular develops single-color genetically encoded biosensors for monitoring signal transduction pathways in living cells. Our sensors can be combined with one another for the simultaneous detection of multiple analytes. Viral packaging means that signaling can be measured in any cell type.

## GPCR Assays in Living Cells

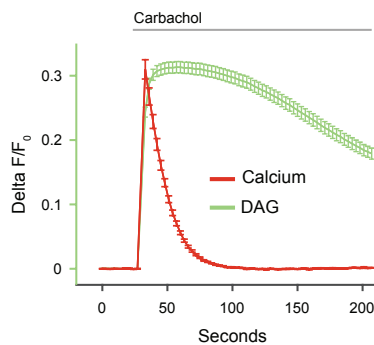
- Gi, Gq, Gs and arrestin pathways.
- Viral delivery to any cell type.
- Easy kinetic analysis for partial and biased agonism.
- No need to add enzyme or co-factors.
- High fluorescence above background.
- Detect on any fluorescence plate reader or imaging system.



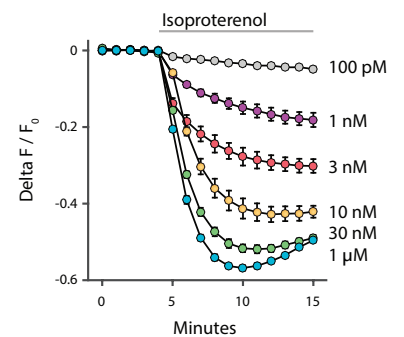
Arrestin – mu opioid receptor



Gi cADDis assay for cAMP

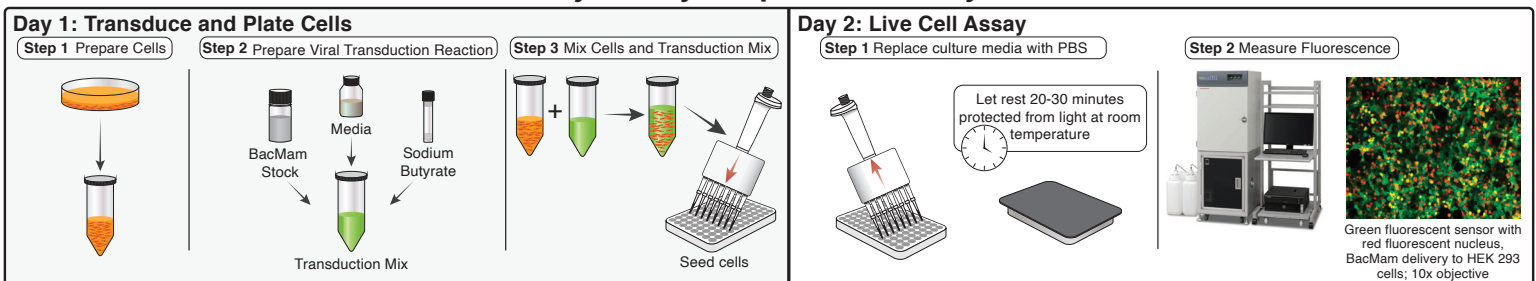


M1 muscarinic receptor - Gq coupled



cADDis assay for cAMP  
β2 adrenergic receptor - Gs coupled

## Easy assay steps - No cell lysis



## Selected References:

Hoare, et al. A kinetic method for measuring agonist efficacy and ligand bias using high resolution biosensors and a kinetic data analysis framework. *Nature Scientific Reports*, Feb 2020.

Hilgendorf, et al. Omega-3 Fatty Acids Activate Ciliary FFAR4 to Control Adipogenesis Cell, November, 2019.

Tewson, et al. Assay for Detecting Gai-Mediated Decreases in cAMP in Living Cells. *SLAS Discovery*. July 10, 2018.

Wray, et al. NMDAR-independent, cAMP-dependent antidepressant actions of ketamine. *Molecular Psychiatry*, 2018.

Almaça, et al. Human beta cells produce and release serotonin to inhibit glucagon secretion from alpha cells. *Cell Reports*, 2016.

Shiget, et al. GLP-1 stimulates insulin secretion by PKC-dependent TRPM4 and TRPM5 activation. *J Clin Invest*. 2015.

Does PKC activation increase the homologous desensitization of μ opioid receptors? *Br J Pharmacy*. 2015.

PKC-dependent Phosphorylation of the H1 Histamine Receptor Modulates TRPC6 Activity. *Cells*. 2014.

Tewson, et al. Simultaneous Detection of Ca<sup>2+</sup> and Diacylglycerol Signaling in Living Cells. *PLoS ONE*. 2012.