



Montana Molecular

Measuring dynamic and real-time cAMP levels using cADDIs, a live-cell indicator for Gs and Gi signaling

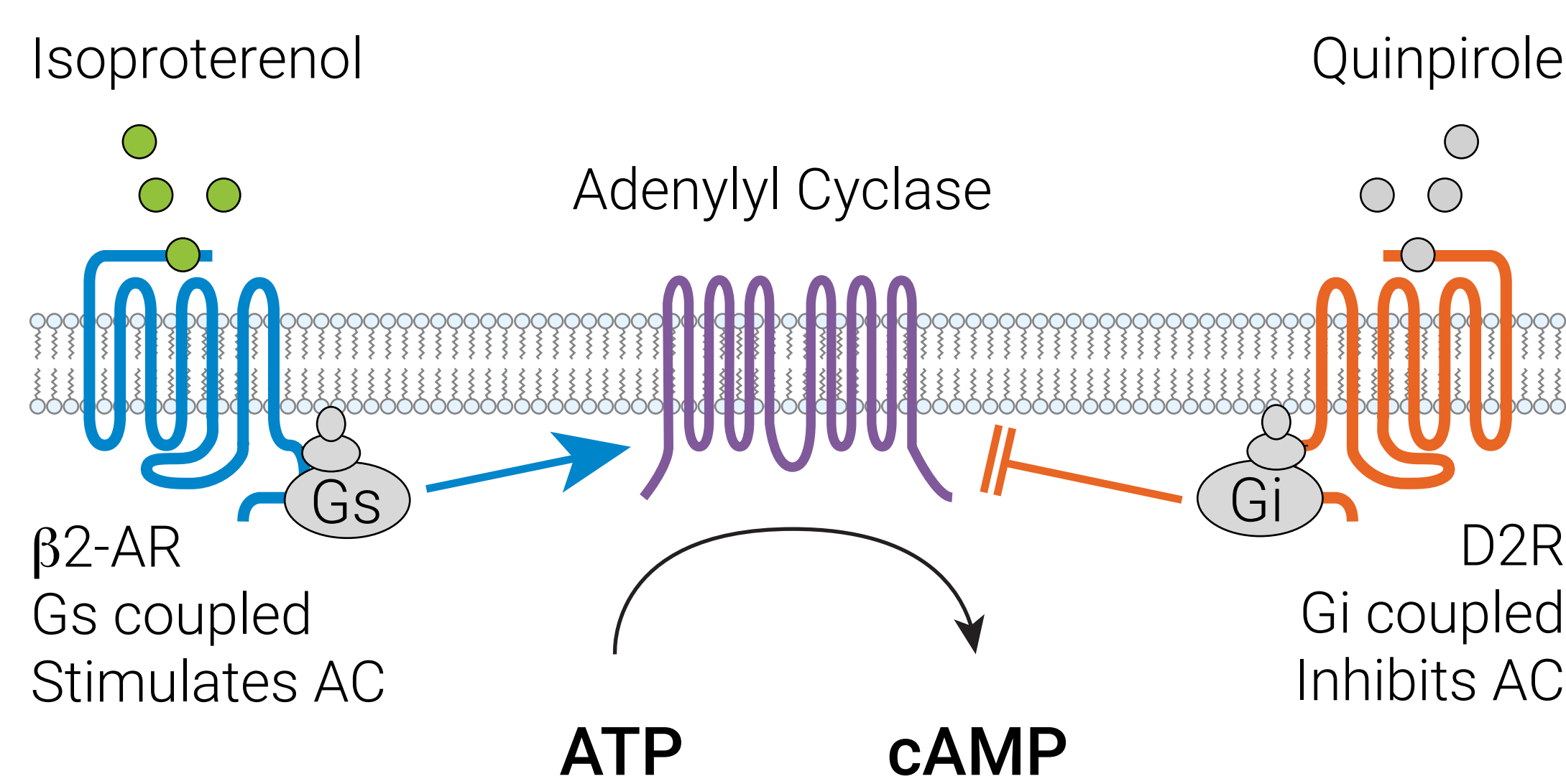
Shane Tillo, Paul Tewson, Scott Martinka, Catherine Berlot, Thom Hughes & Anne Marie Quinn



MM cADDIs Sensors for Drug Discovery

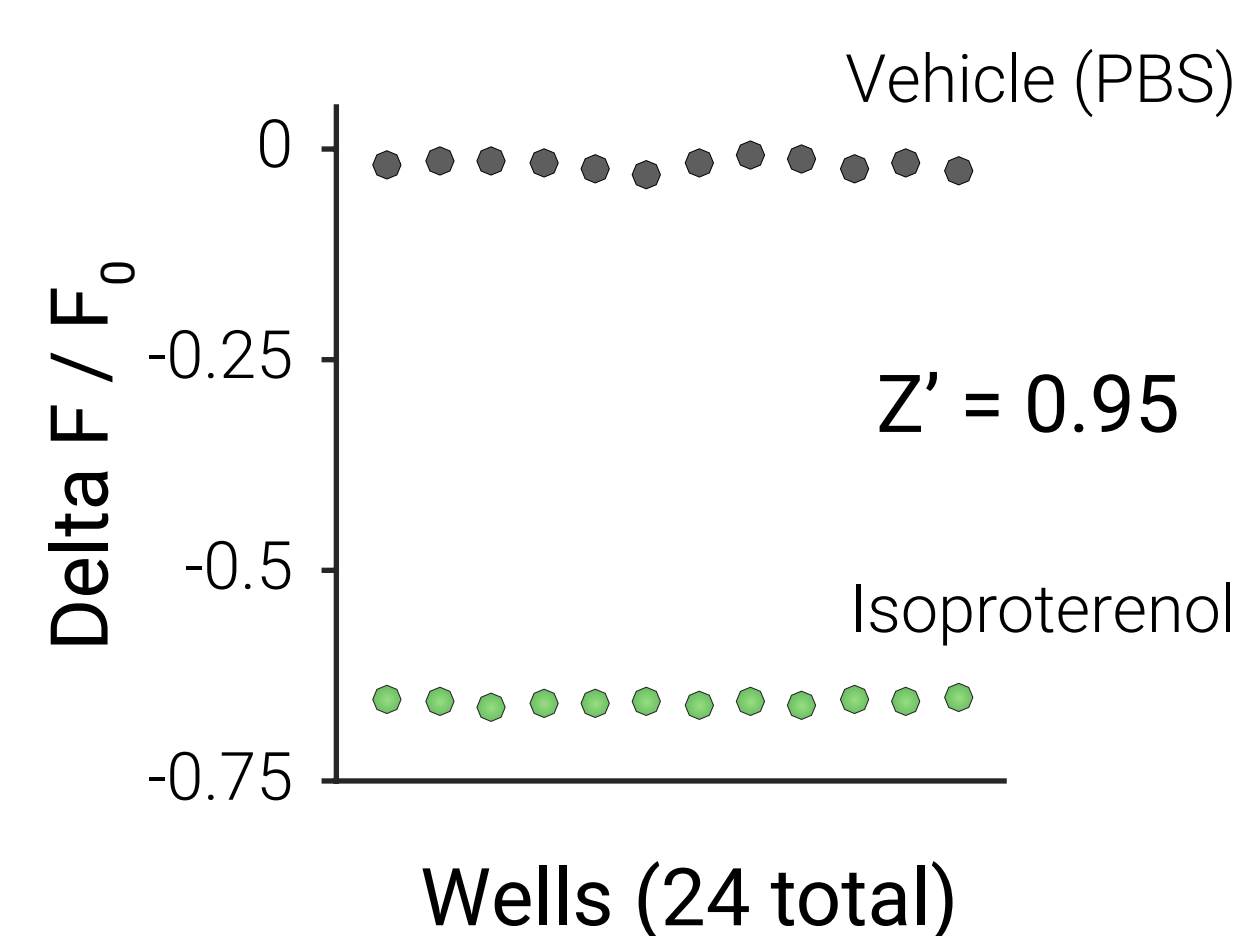
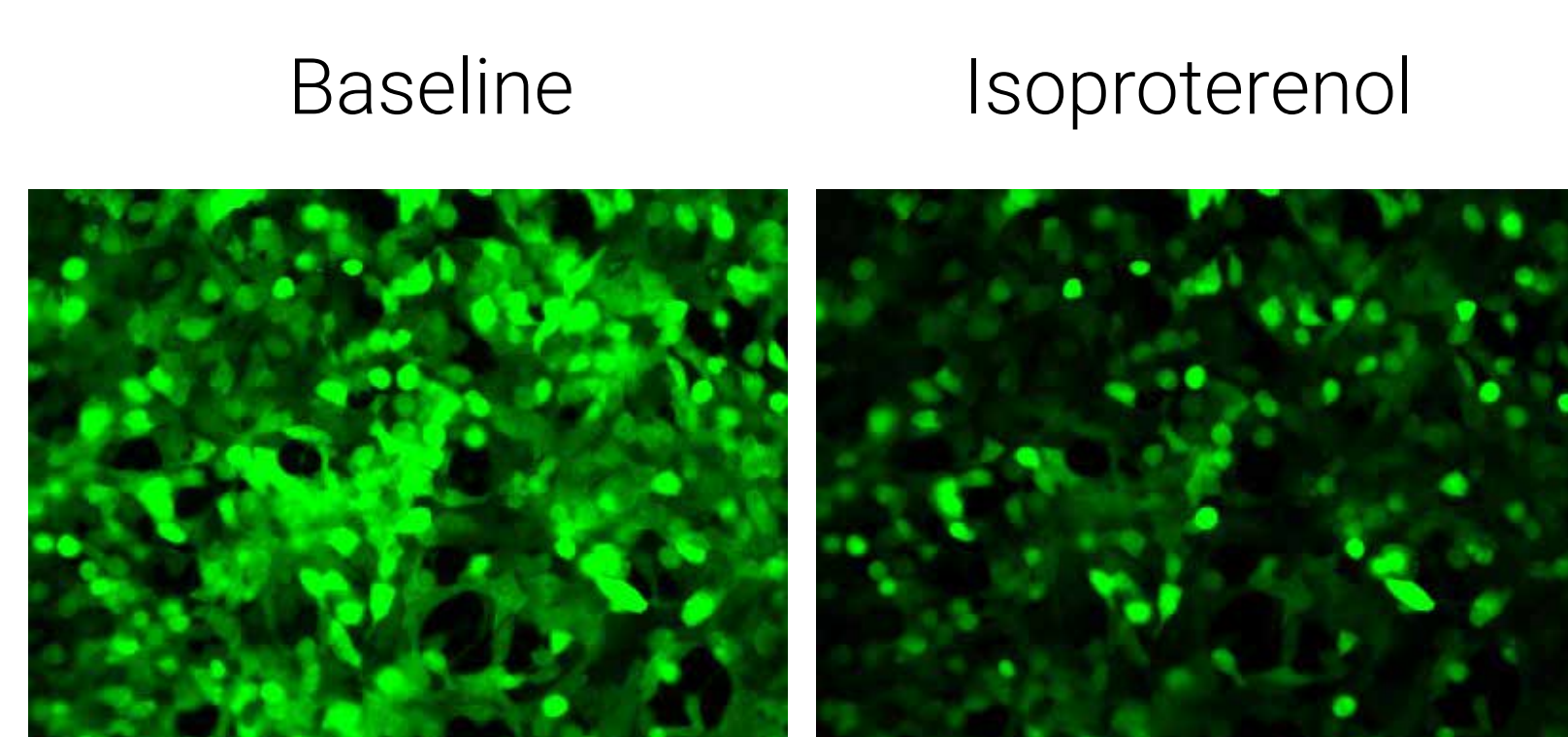
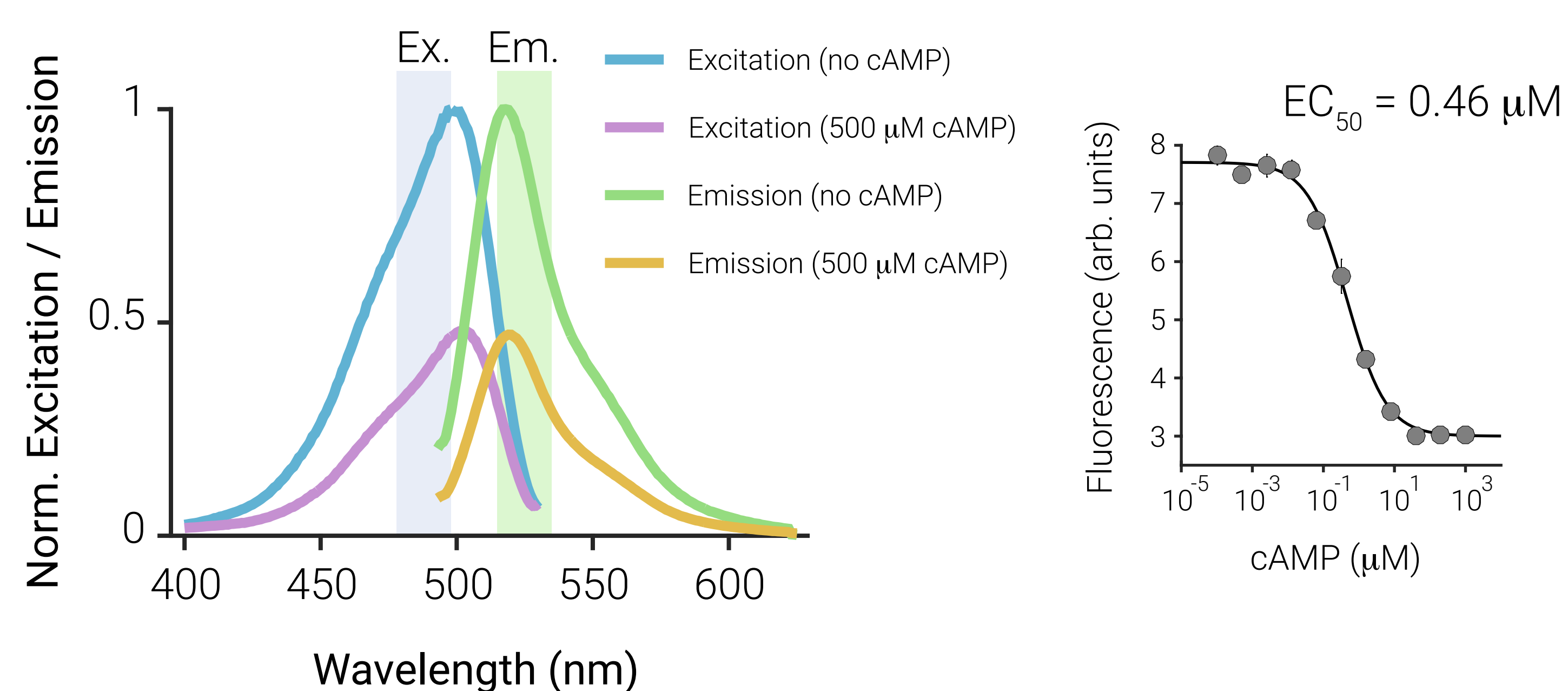
Real-Time cAMP Detection

Here, we introduce an improved fluorescent cAMP indicator, cADDIs, capable of detecting dynamic changes in cAMP concentration in living cells. We have applied cADDIs to use in a novel Gi assay that, for the first time, directly reports Gi mediated reductions in cAMP concentration, across a variety of receptors, on standard automated fluorescence plate readers.



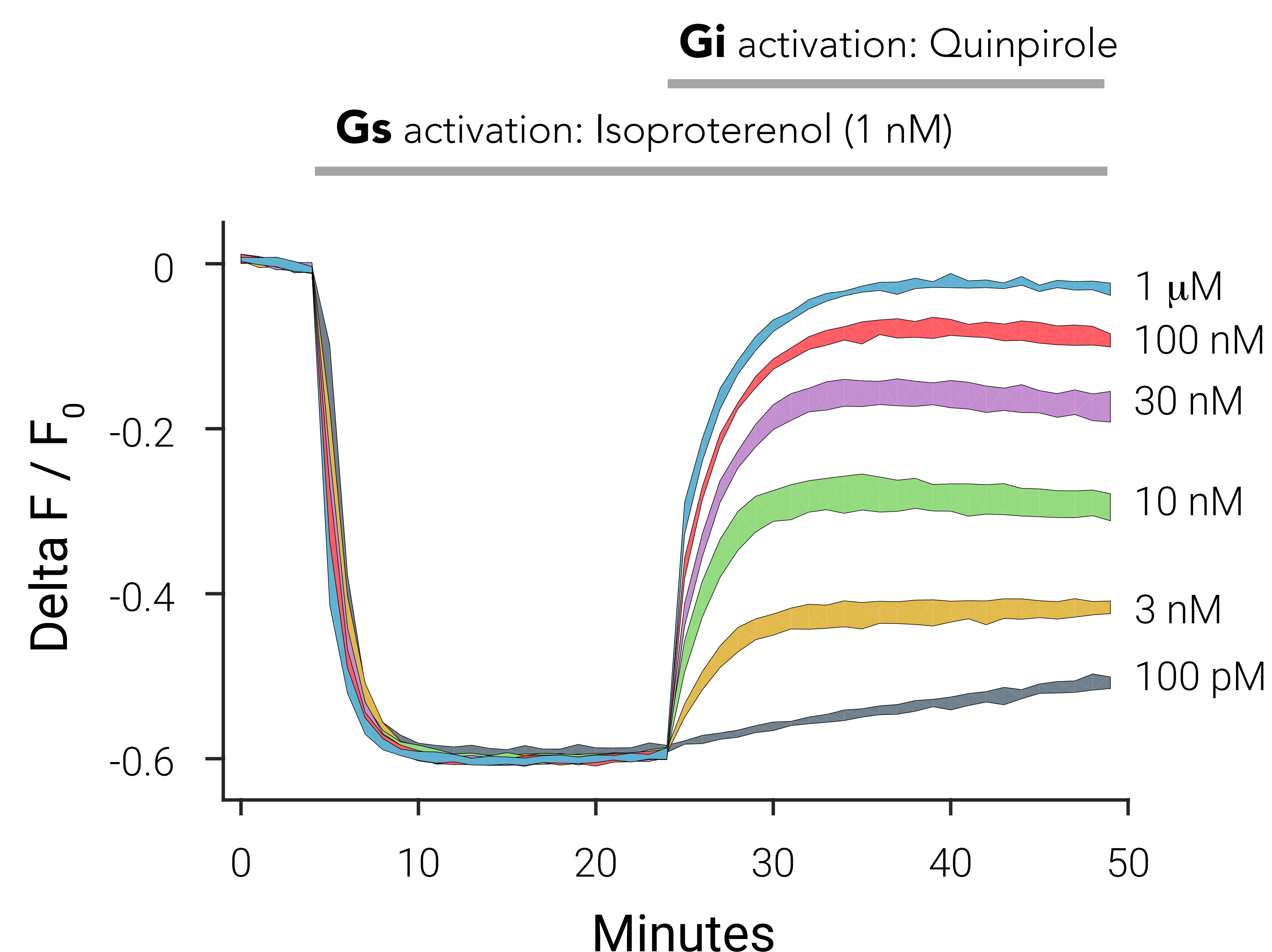
cADDIs cAMP Indicator

- Genetically encoded, single fluorescence emission
- Packaged in a variety of viral vectors for use in any cell type
- Decrease in fluorescence signal upon binding cAMP
- Easy to use on standard fluorescence microscopes and automated fluorescence plate readers



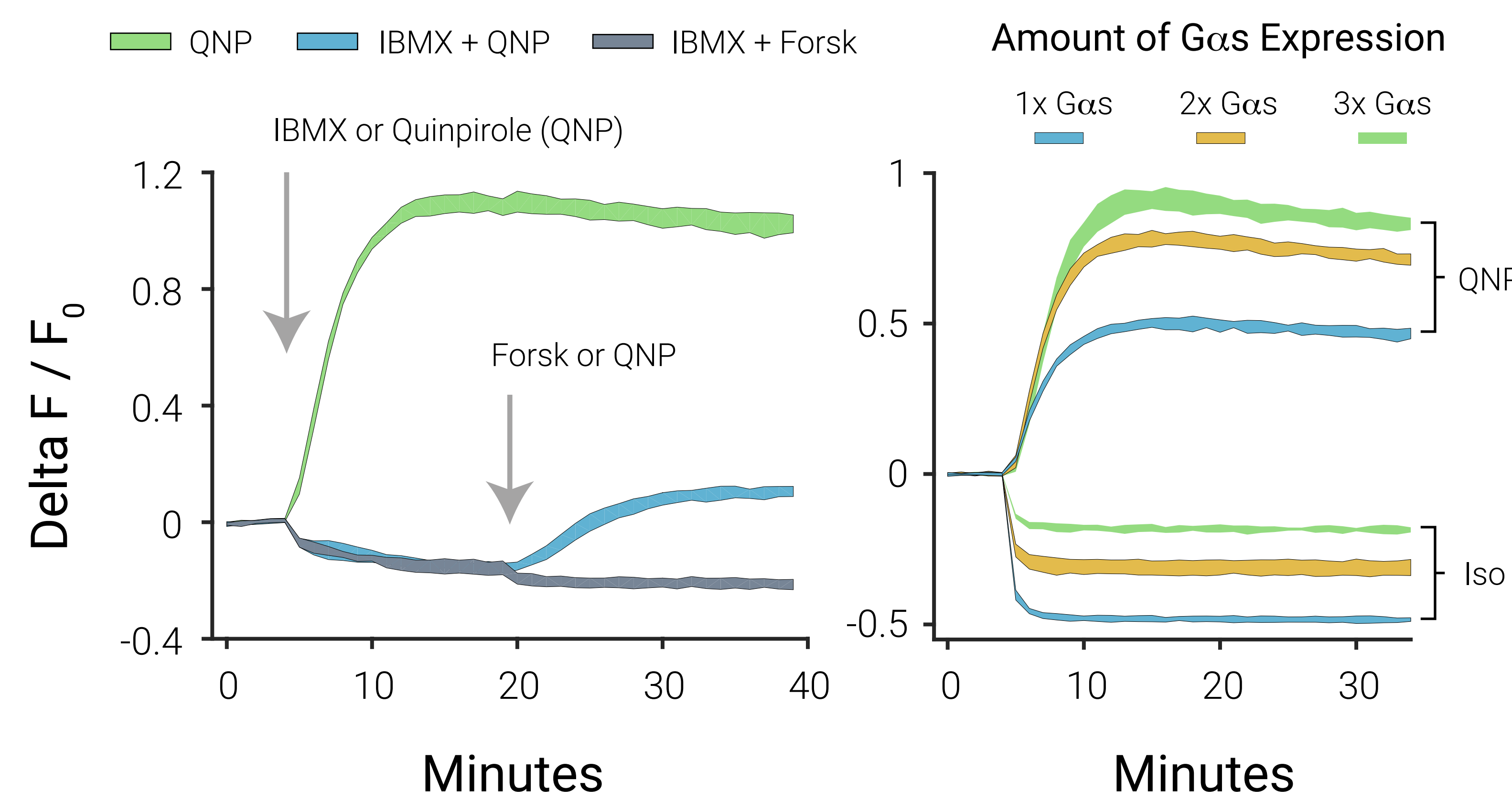
A New Way to Detect Gi Signaling

- Simple, No cell lysis, No FRET, Fully reversible
- Real-time detection of dynamic cAMP levels
- All data obtained on Biotek Synergy MX plate reader

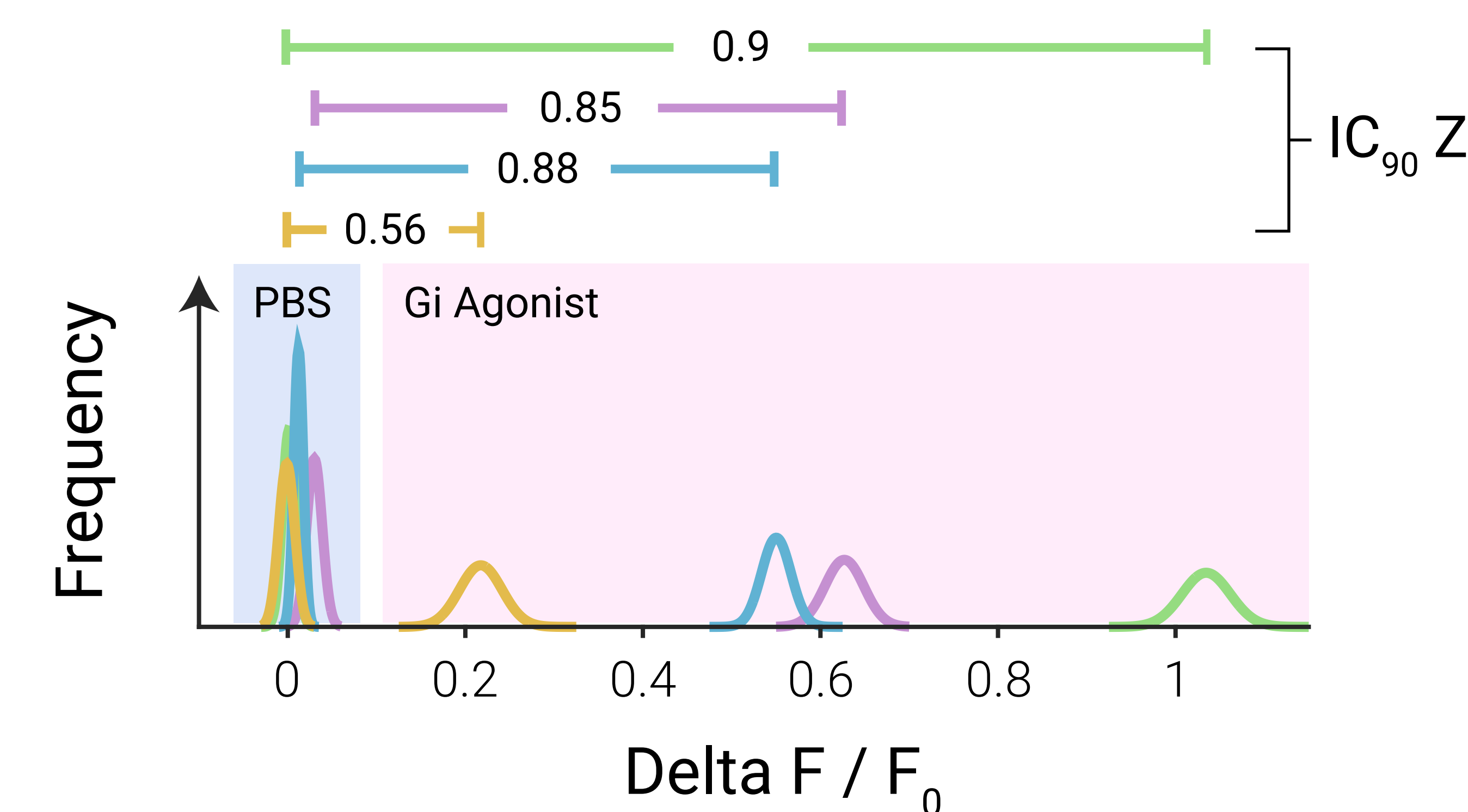
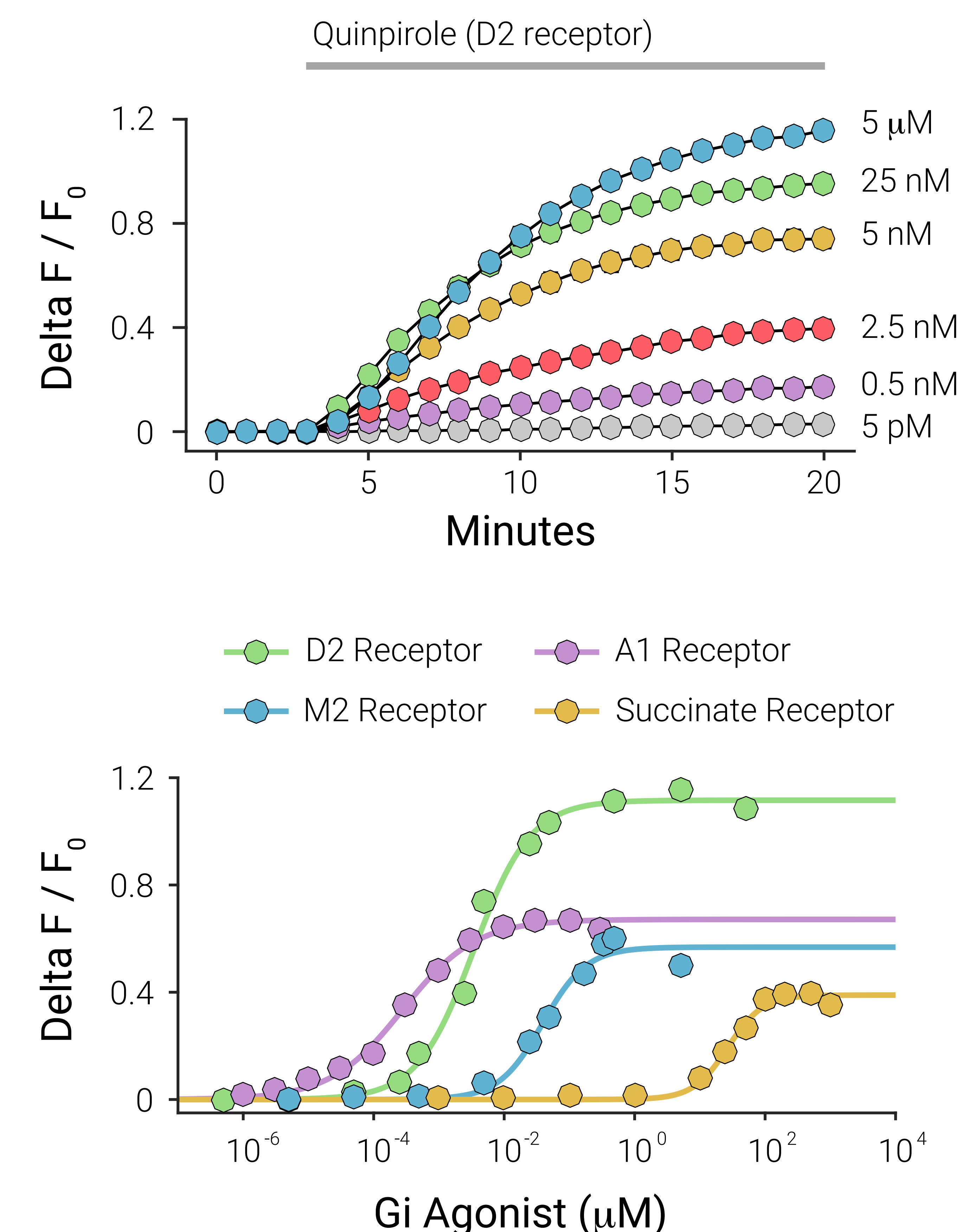


First ever direct readout of Gi mediated reductions in cAMP

- Tuned expression of constitutively active Gαs raises basal levels of cAMP, without saturating cADDIs2
- No pre-treatment with Gs agonists, Forskolin, or IBMX
- Reveal the activity of Phosphodiesterases



Direct, Real-Time Gi Detection



Acknowledgements:

Montana Molecular sensors made possible by:
 NSF SBIR IIP-1430878
 NIH SBIR R44 NS082222
 Montana Dept. of Commerce MBRCT: #15-22