

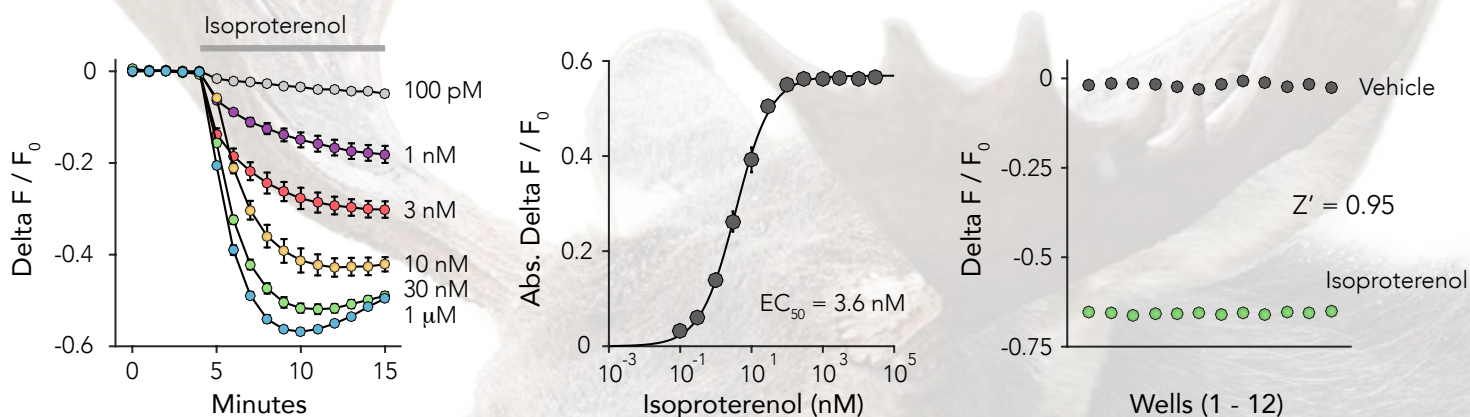
Introducing cADDIs: The simplest assay in the world for detecting cAMP

Measure cAMP signaling in real time, in the living cells of your choice.

Save time and money by switching to cADDIs, a simple and robust assay for Gs and Gi signaling, without the need for co-factors, cell lysis, or complex liquid handling. cADDIs is genetically-encoded and packaged in BacMam for easy transduction into a wide variety of cell types, including primary and iPSC-derived cell lines. Simply add cADDIs to your cells and measure robust signals the very next day on standard fluorescence microscopes and automated plate readers.

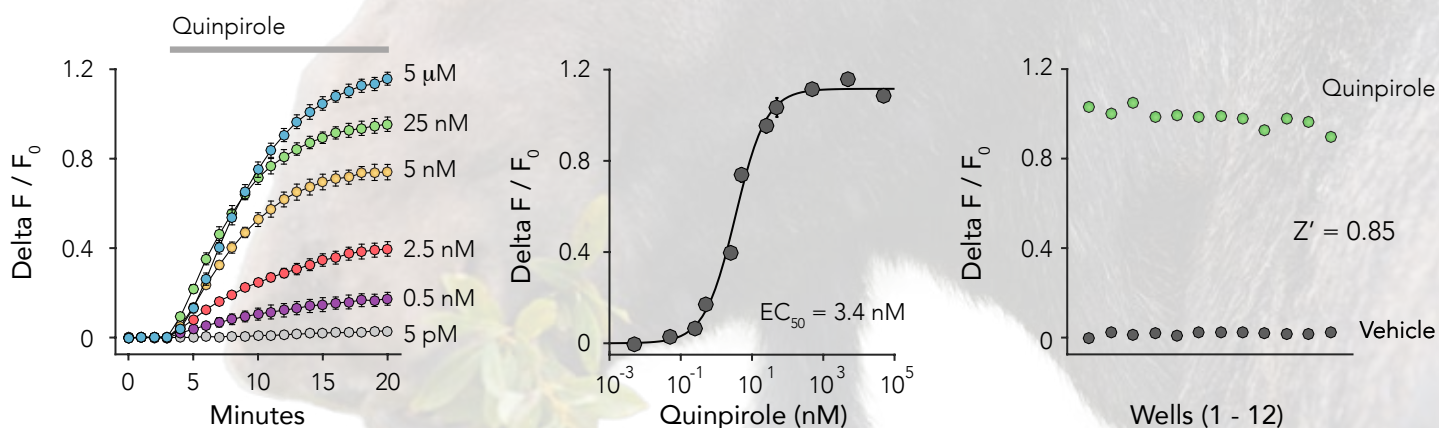
Gs detection: Kinetics. Dose responses. Bright and reliable signals.

Measure kinetics and dose responses of Gs-mediated cAMP production with large and consistent fluorescence changes. cADDIs is extremely bright with remarkable signal to noise ($Z' > 0.95$). Below: cADDIs in live HEK293 cells stimulated with isoproterenol, a selective agonist of β -adrenergic receptors.



Gi detection: Forskolin-free

cADDIs detects Gi-mediated decreases in cAMP production, in real time, without the need for forskolin or IBMX. The cADDIs Gi assay yields Z' values > 0.85 . Forego the forskolin and measure biologically relevant Gi signaling using cADDIs. Below: cADDIs in live HEK293 cells stimulated with quinpirole, a selective agonist of dopaminergic D2 receptors.



Discover cADDIs today

Contact: info@montanamolecular.com