

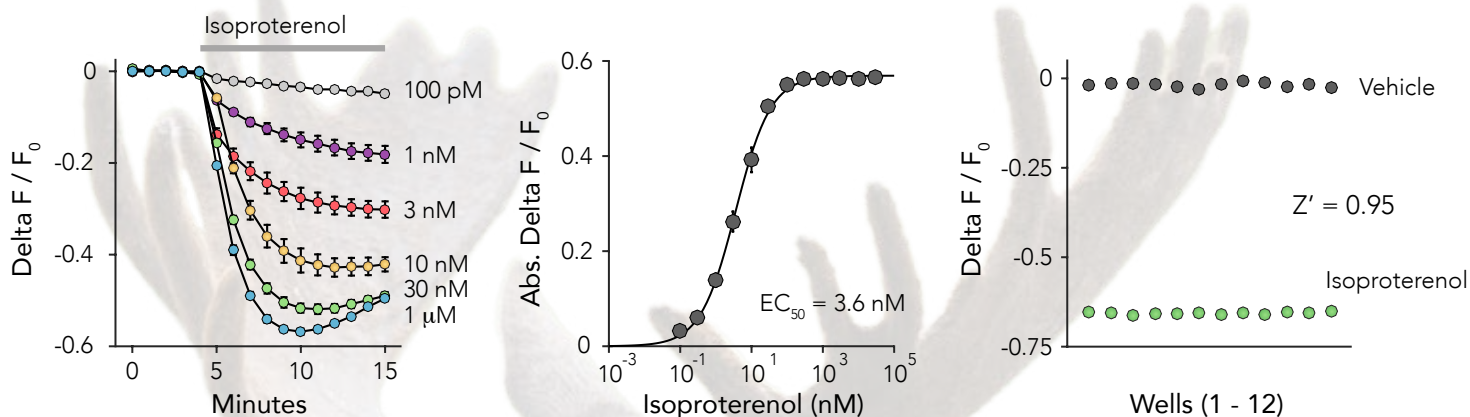
# Introducing cADDis: The simplest assay in the world for detecting cAMP

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.

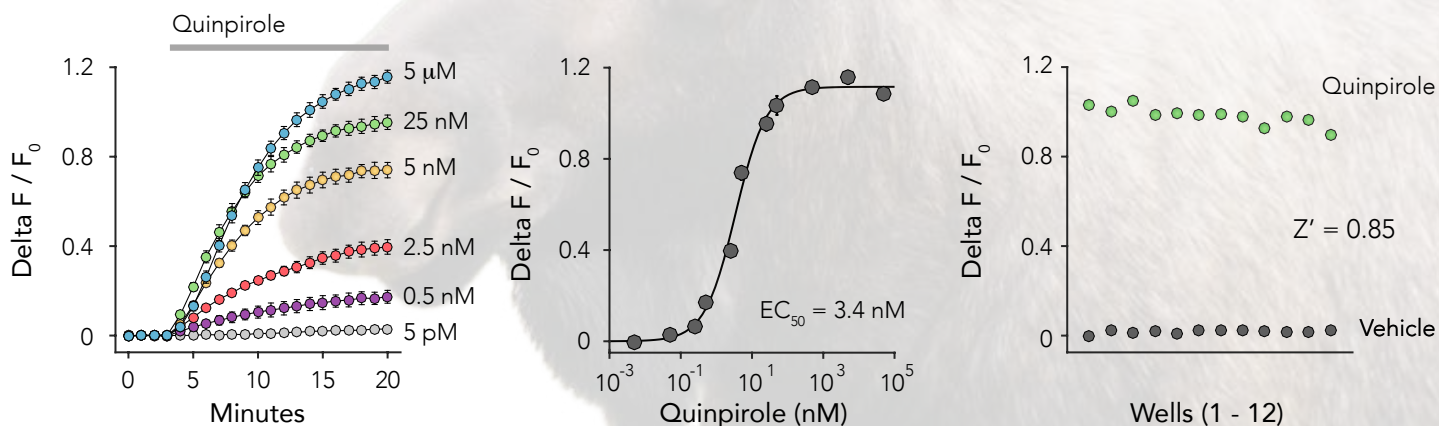
## Gs detection: Kinetics. Dose responses. Bright and reliable signals.

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  $\beta$ -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](mailto:info@montanamolecular.com)