

Monitoring the Activity of GPCR Modulated by Lipid or Free Fatty Acid Agonists

ABSTRACT

We have recently demonstrated that the bioluminescent cAMP-Glo™ Assay can track direct activation of adenylate cyclase by forskolin as well as agonists and antagonists that modulate GPCRs by monitoring the level of intracellular cAMP in the cell with receptors that modulate Gas (Dopamine D1) and those that modulate Gai (Dopamine D2). Here we extend the application of the cAMP-Glo™ Assay using receptors that are modulated by lipid and free fatty acid agonists. These receptors are of interest since little is known about them and there is conflicting information in the available literature.

Cell Notes 23, 13–16.

Said Goueli and Kevin Hsaio

Promega Corporation

Publication Date: 2009



Download Article

(249 KB)

HOW TO CITE THIS ARTICLE

Scientific Style and Format, 7th edition, 2006

Goueli, S. and Hsaio, K. Monitoring the Activity of G Protein-Coupled Receptors (GPCR) Modulated by Lipid or Free Fatty Acid Agonists. [Internet] 2009. [cited: year, month, date]. Available from: <http://www.promega.com/resources/pubhub/cellnotes/monitoring-the-activity-of-gpcrs--modulated-by-lipids-or-free-fatty-acids/>

American Medical Association, Manual of Style, 10th edition, 2007

Goueli, S. and Hsaio, K. Monitoring the Activity of G Protein-Coupled Receptors (GPCR) Modulated by Lipid or Free Fatty Acid Agonists. Promega Corporation Web site. <http://www.promega.com/resources/pubhub/cellnotes/monitoring-the-activity-of-gpcrs--modulated-by-lipids-or-free-fatty-acids/> Updated 2009. Accessed Month Day, Year.