

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT NPFF2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CG1364

Lot Number: CG1364-101215

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gaqi5

Transfection: Expression vector containing full-length human NPFF1 cDNA (GenBank Accession Number NM_004885.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: Neuropeptide FF receptor 2 (NPFF2) belongs to the rhodopsin family of G protein-coupled receptors and found across the central nervous system. It is believed to regulate endogenous opioid systems in the brain and spinal cord. The neuropeptide is considered a prospective target to decrease negative effects of chronic opioid use.

Application: Functional assays

Figure 1

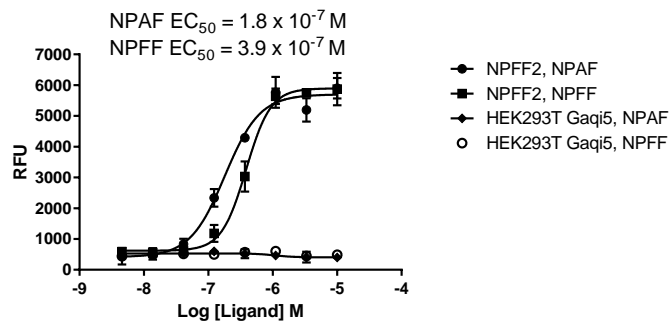


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Bonini, James et al. "Identification and Characterization of Two G Protein-coupled Receptors for Neuropeptide FF." *The Journal of Biological Chemistry*. Vol. 275, No. 50, pp. 39324-39331. Dec 2000

Elshourbagy, Nabil et al. "Receptor for the Pain Modulatory Neuropeptides FF and AF is an Orphan G Protein-coupled Receptor." *The Journal of Biological Chemistry*. Vol. 275, No. 34, pp. 25965-25971. Aug 2000

Mollereau, Catherine et al. "Agonist and Antagonist Activities on Human NPFF₂ Receptors of the NPY Ligands GR231118 and BIBP3226." *British Journal of Pharmacology* 133.1 (2001): 1-4. *PMC*. Web. 5 Mar. 2015.

Moulédous, Lionel et al. "Opposite control of body temperature by NPFF1 and NPFF2 receptors in mice." *Neuropeptides* 44.5: 453 - 456.

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.