

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT NPFF1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1363

Lot Number: C1363-061814

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

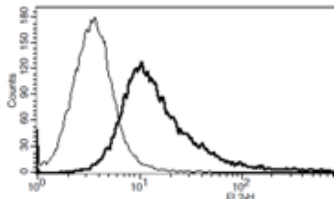
Transfection: Expression vector containing full-length human NPFF1 cDNA (GenBank Accession Number NM_022146.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: In Progress

Figure 2



Background: Neuropeptide FF receptor 1 (NPFF1) acts as a modulator of endogenous opioid function. NPFF1 is found across the central nervous system. It is widely expressed in the hypothalamus and limbic system. Recent studies performed on NPFF1 receptors in mice implicate the receptor as an important regulator of body temperature.

Application: Functional assays

Figure 1

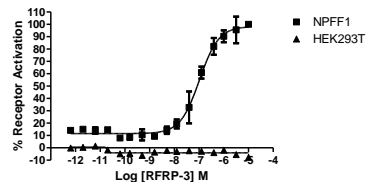


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

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.