

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT NPY2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: CG1274-1

Lot Number: CG1274-1-022008

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gαq5

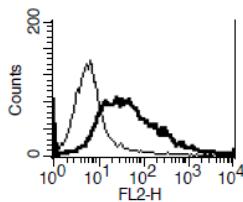
Transfection: Expression vector containing full-length human NPY2R cDNA (GenBank Accession Number NM_00091) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM-F12, 10% FBS, 10 µg/mL puromycin, 150 µg/mL hygromycin

Stability: Stable after two months continuous growth

Figure 3



Background: NPY2R (Y2) belongs to the family of G protein-coupled receptors. It is expressed at high levels in central nerve system, heart, ileum and colon. NPY2R shows high affinity for neuropeptide Y (NPY), peptide YY (PYY), and a fragment of NPY including amino acids 13 to 36. The rank order of affinity of this receptor for pancreatic polypeptides is PYY > NPY > PYY (3-36) > NPY (2-36) > [I¹-31, Gln-34] PP > [Leu-31, Pro-34] NPY > PP, [Pro-34] PYY and NPY free acid. NPY exhibits a variety of potent central and peripheral effects including those on feeding, memory, blood pressure, and intestinal secretions. NPY2R has been shown to contribute to obesity.

Application: Functional assays

Figure 1

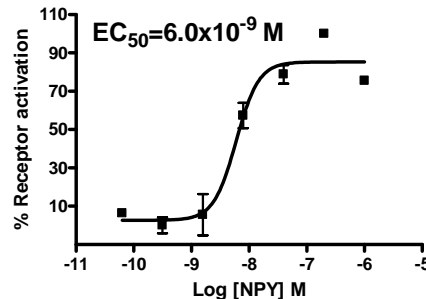
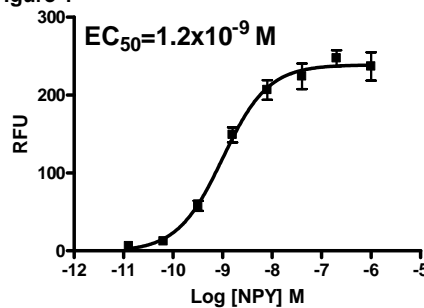


Figure 2

Figure 1. Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent cAMP accumulation upon treatment with ligand. **Figure 3.** 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:

Gerald *et al.* (1995) Expression cloning and pharmacological characterization of a human hippocampal neuropeptide Y/peptide YY Y2 receptor subtype. *J Biol Chem* 270:26758-26761.

Lutz *et al.* (1997) Neuropeptide Y receptor genes on human chromosome 4q31-q32 map to conserved linkage groups on mouse chromosomes 3 and 8. *Genomics* 41:498-500.

Rose *et al.* (1995) Cloning and functional expression of a cDNA encoding a human type 2 neuropeptide Y receptor. *J Biol Chem* 270:22661-22664.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.