

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT NPY4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DCG1300-1

Lot Number: DCG1300-1-060122

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1 Gaq15

Transfection: Expression vector containing full-length human Y4 cDNA (GenBank Accession Number AY268432) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM-F12, 10% FBS

Stability: Stable for 1 – 2 days after thawing

Background: NPY4R (Y4) has been reported to be expressed at highest levels in brain, coronary artery and ileum, low levels in pancreas and kidney. It is also been detected in colon and small intestine and isolated from human colon and head/neck libraries. It is one of the receptors for neuropeptide Y and peptide YY. The rank order of affinity of this receptor for pancreatic polypeptides is PP, PP (2-36) and [Ile-31, Gln-34] PP > [Pro-34] PYY > PYY and [Leu-31, Pro-34] NPY > NPY > PYY (3-36) and NPY (2-36) > PP (13-36) > PP (31-36) > NPY free acid. Ligand: Neuropeptide Y, Peptide YY.

Application: Functional assay

Figure 1

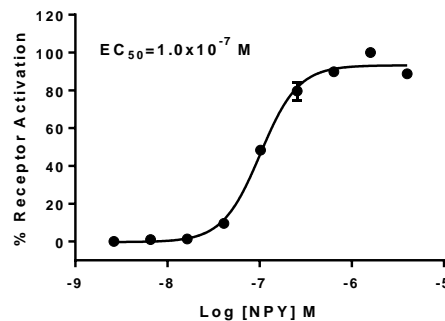


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Bard et al. (1995) Cloning and functional expression of a human Y4 subtype receptor for pancreatic polypeptide, neuropeptide Y, and peptide YY. *J Biol Chem* 270:26762-26765.

Yan et al. (1996) Cloning and functional expression of cDNAs encoding human and rat pancreatic polypeptide receptors. *Proc Natl Acad Sci USA* 93:4661-4665.

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.