

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT Y5 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1275a

Lot Number: C1275a-082510

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length Human NPY5R cDNA (GenBank Accession NM_006174) with FLAG-tag sequence at the 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: NPY5R encodes 456-amino acid protein with 7 putative transmembrane domains. It has shown to be expressed in the brain and hypothalamus. Y5 receptors' pharmacological profile correlates it with peptides known to elicit a feeding response. It has been suggested to have effects on psychomotor activity, food intake, regulation of central endocrine secretion, and potent vasoactive effects on the cardiovascular system.

Application: Functional assays

Figure 1

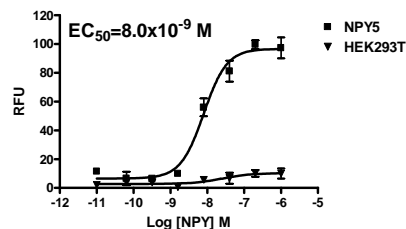


Figure 2

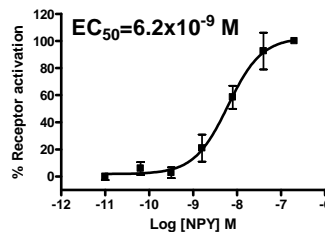


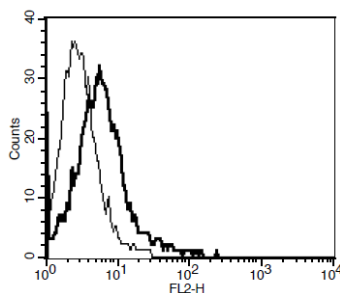
Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 3.** Receptor expression 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.* (1996) A receptor subtype involved in neuropeptide-Y induced food intake. *Nature* 382:168-171.

Hu *et al.* (1996) Identification of a novel hypothalamic neuropeptide Y receptor associated with feeding behavior. *J Biol Chem* 271:26315-26319.

Figure 3



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.