

MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT NPSR1 RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1355-1

Lot Number: C1355-1-021909

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length mouse Npsr1 cDNA (GenBank Accession Number NM_175678)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 100 μ g/mL hygromycin, 100 μ g/mL G418

Stability: Stable after minimum of two months continuous growth

Data sheet

Background: Neuropeptide S (NPS) receptor 1 is also known as GPR154 or G protein-coupled receptor for asthma susceptibility (GPRA). NPS receptor is widely expressed in the brain, with highest levels found in hypothalamus, amygdala, endopiriform nucleus, cortex, subiculum and nuclei of the thalamic midline. Central administration of NPS promotes behavioral arousal and suppresses all stages of sleep in rodents. Furthermore, NPS was found to produce anxiolytic-like effects in a battery of four different tests that measure behavioral responses of rodents to novelty or stress.

Application: Functional assays

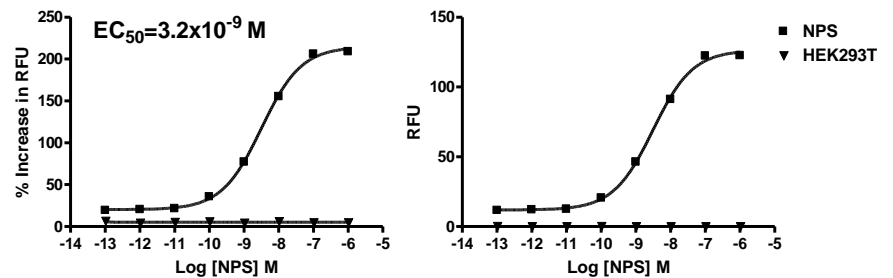


Figure legend: Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Xu *et al.* (2004) Neuropeptide S: a neuropeptide promoting arousal and anxiolytic-like effects. *Neuron* 43:487-497.

Laitinen *et al.* (2004) Characterization of a common susceptibility locus for asthma-related traits. *Science* 304:300-304.

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. June 2005

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