

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT NMUR1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1122

Lot Number: C1122-091505

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 NMUR1 cDNA (GenBank Accession Number NM_006056) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum of two months continuous growth

Background: NMUR1, or FM-3 or GPR66, is a receptor for the neuromedin U, a neuropeptide that has been implicated in physiological roles, including the regulation of feeding, anxiety, pain, blood flow, and smooth muscle contraction. The NMUR1 is abundantly expressed in peripheral tissues such as pancreas, testis and small intestine.

Application: Functional assays

Figure 1

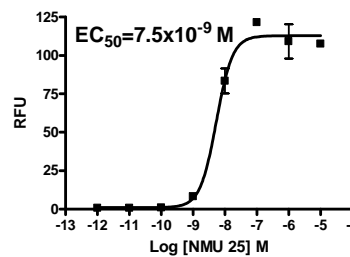


Figure 2

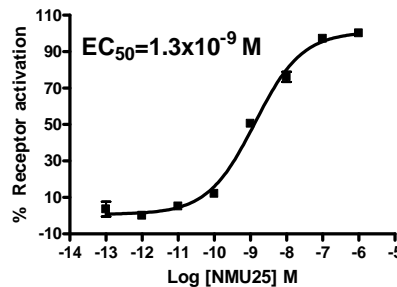


Figure 3

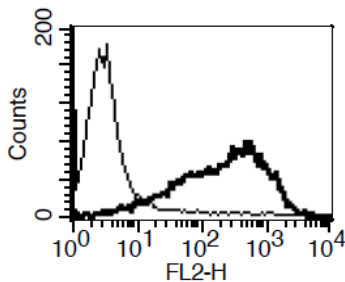


Figure 1. Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent accumulation of intracellular IP1 upon treatment with ligand, measured with IP-one Tb kit (Cisbio 62IPAPEC). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

- Howard *et al.* (2000) Identification of receptors for neuromedin U and its role in feeding. *Nature* 406:70-74.
- Kojima *et al.* (2000) Purification and identification of neuromedin U as an endogenous ligand for an orphan receptor GPR66 (FM3). *Biochem Biophys Res Commun* 276:435-438.

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.