

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

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1122

**Lot Number:** C1122-091505

**Quantity:** 1 vial (2 x 10<sup>6</sup>) 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

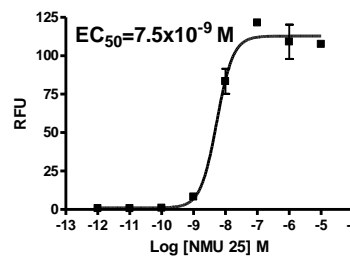
**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin

**Stability:** Stable after minimum of two months continuous growth

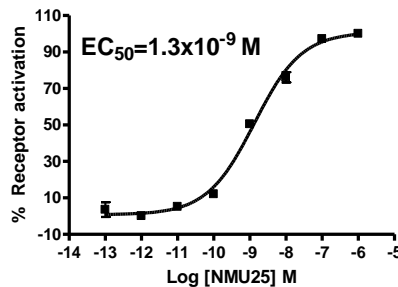
**Background:** NMUR1, or FM-3 or GPR66, is a receptor for the neuropeptide U, a neuropeptide that has been implicated in physiological roles, including the regulation of feeding, anxiety, pain, blood flow, and smooth muscle contraction. The NMUR1 is abundantly expressed in peripheral tissues such as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

**Application:** Functional assays

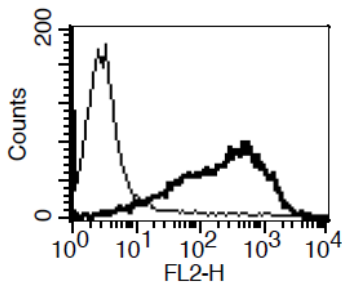
**Figure 1**



**Figure 2**



**Figure 3**



**Figure 1.** Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Dose-dependent accumulation of intracellular IP1 upon treatment with ligand, measured with IP-one Tb kit. **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 neuropeptide U and its role in feeding. *Nature* 406:70-74.
- Kojima *et al.* (2000) Purification and identification of neuropeptide U as an endogenous ligand for an orphan receptor GPR66 (FM3). *Biochem Biophys Res Commun* 276:435-438.

**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.