

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT NMUR1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1122

Lot Number: DC1122-041619

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2

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

Stability: 1-2 days after thawing

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

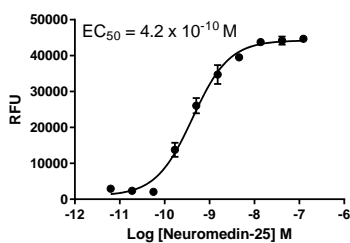


Figure 1. Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.