

## MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT NMU2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DC1213

**Lot Number:** DC1213-041619

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

**Freeze Medium:** Cellbanker 2

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human NMUR2 cDNA (one sense mutation, which has been noted in literature, in comparison to GenBank Accession Number NM\_020167.3) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

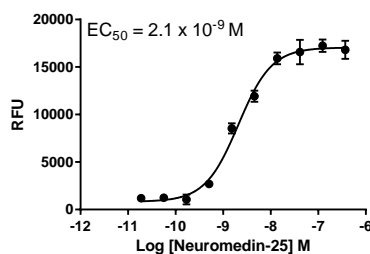
**Propagation Medium:** DMEM, 10% FBS

**Stability:** 1-2 days after thawing

**Background:** NMUR2 is the second receptor for the neuromedin U, which is a neuropeptide and has been implicated in physiological roles, including the regulation of feeding, anxiety, pain, blood flow and smooth muscle contraction. NMUR2 is expressed predominantly in the central nervous system. Central administration of neuromedin U (NMU) suppresses food intake acting through the NMU-2 receptor (NMU2R), which is expressed in the hypothalamus.

**Application:** Functional assays

**Figure 1**



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

#### References:

Bhattacharyya *et al.* (2004) Studies of the neuromedin U-2 receptor gene in human obesity: evidence for the existence of two ancestral forms of the receptor. *J Endocrinol* 183:115-120.

Brighton *et al.* (2004) Signaling and ligand binding by recombinant neuromedin U receptors: evidence for dual coupling to Galphaq/11 and Galphai and an irreversible ligand-receptor interaction. *Mol Pharmacol* 66:1544-1556.

**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.