

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT MRGPRX2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1257

**Lot Number:** C1257-073005

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T Gaq15

**Transfection:** Full-length human MRGPRX2 cDNA (GenBank Accession Number NM\_054030)

**Recommended Storage:** Liquid nitrogen upon receiving

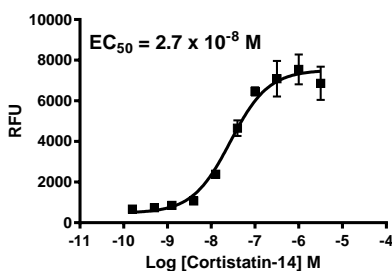
**Propagation Medium:** DMEM, 10% FBS, 200  $\mu$ g/mL hygromycin, 0.625  $\mu$ g/mL puromycin

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

**Background:** MRGPRX2 (MAS-related GPR member X2) is a receptor for cortistatin. It is probably involved in the function of nociceptive neurons and regulation of nociceptor function and/or development, including the sensation or modulation of pain. Cortistatin-14, a high potency agonist at the receptor, has biological functions including sleep regulation, locomotor activity and cortical function.

**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:

Robas *et al.* (2003) MrgX2 is a high potency cortistatin receptor expressed in dorsal root ganglion. *J Biol Chem* 278:44400-44404.

Kamohara *et al.* (2005) Identification of MrgX2 as a human G-protein-coupled receptor for proadrenomedullin N-terminal peptides. *Biochem Biophys Res Commun* 330:1146-1152.

**FOR RESEARCH USE ONLY.**

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