

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT MRGX1/SNSR4 RECEPTOR**

**PRODUCT INFORMATION**

**Catalog Number:** C1256

**Lot Number:** C1256-061505

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

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

**Host cell:** CHO dhfr<sup>-</sup> Gaqi5

**Transfection:** Full-length human MRGX1 cDNA (GenBank Accession Number NM\_147199)

**Recommended Storage:** Liquid nitrogen upon receiving

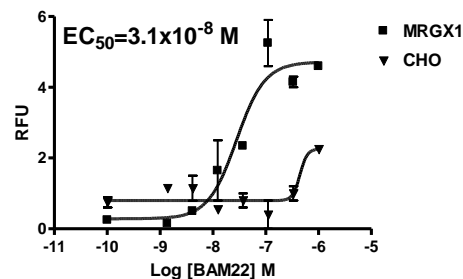
**Propagation Medium:** Alpha-MEM, 10% FBS, 100 µg/mL hygromycin, 800 µg/mL G418

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

**Data sheet**

**Background:** MRGX1 (MAS-related GPR member X1) is also known as SNSR4 (sensory neuron-specific G-protein-coupled receptor 4). It can be potently activated by enkephalins including BAM22 and BAM(8-22). MRGX1 receptor is expressed solely in small diameter primary sensory neurons. This restricted expression pattern is of considerable therapeutic interest because small nociceptors transmit chronic pain messages.

**Application:** Functional assays



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

**References:**

Chen and Ikeda (2004) Modulation of ion channels and synaptic transmission by a human sensory neuron-specific G-protein-coupled receptor, SNSR4/mrgX1, heterologously expressed in cultured rat neurons. *J Neurosci* 24:5044-5053.

Ahmad and Dray (2004) Novel G protein-coupled receptors as pain targets. *Curr Opin Investig Drugs* 5:67-70.

**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.