

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT MRGX2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1257

Lot Number: C1257-073005

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gaqi5

Transfection: Full-length human MRGX2 cDNA (GenBank Accession Number NM_054030)

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum of two months continuous growth

Background: MRGX2 (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

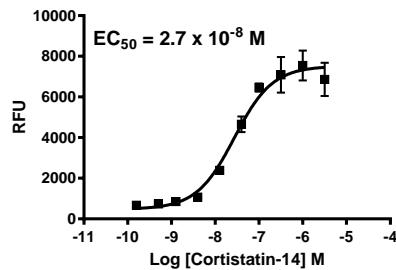


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.