

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT MRGPRX2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1257b

Lot Number: MC1257b-101624

Quantity: Available 1, 5, or 10mg per vial

Protein concentration: 15.12 mg/mL

Packaging Buffer: 20mM Gly-Gly, 1 mM MgCl₂, 25mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Full-length human MRGPRX2 cDNA (GenBank Accession Number NM_054030) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Background: MRGPRX2 (MAS-related GPR member X2) is widely known as 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. Recent studies have found that MrgX2 is also a potential human PAMP-12 receptor that acts by inhibiting of the forskolin-elevated cAMP accumulation and regulates catecholamine secretion from adrenal glands.

Application: Functional assays

Figure 1

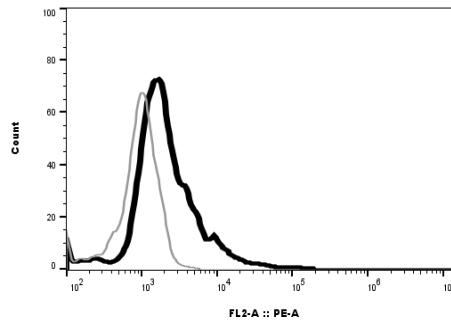


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.