

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT S1P4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MCG1052-1

Lot Number: MCG1052-1-041321

Quantity: 1 vial (4.6 mg/mL),
1mg

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

Host cell: CHO-K1 Gaqi5

Transfection: Expression vector
containing full-length human S1P4
cDNA (GenBank Accession Number
NM_003775) with FLAG tag sequence
at N-terminus

Recommended Storage: Liquid
nitrogen upon receiving. Avoid
repeated freeze-thaw.

Background: S1P4 receptor or Endothelial Differentiation Gene-6 (EDG-6) is receptor for sphingosine-1-phosphate (S1P). S1P is a bioactive lipid produced from the metabolism of sphingomyelin. It is an important constituent of serum and regulates cell growth, survival, migration, differentiation and gene expression. Unlike other members of S1P receptor, which are widely expressed, S1P4 exhibits lymphoid tissue-specific expression. EDG-6 has been implicated in regulation of cell shape and motility.

Figure 1

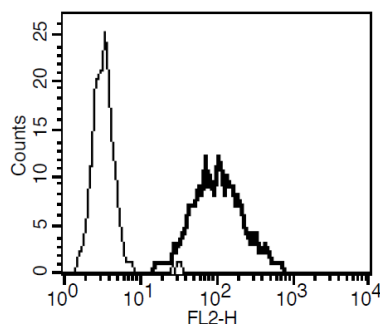


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody.

References:

Clemens *et al.* (2004) Synthesis of benzimidazole based analogues of sphingosine-1-phosphate: discovery of potent, subtype-selective S1P4 receptor agonists. *Bioorg Med Chem Lett* 14:4903-4906.

Takuwa *et al.* (2001) Subtype-specific, differential activities of the EDG family receptors for sphingosine-1-phosphate, a novel lysophospholipid mediator. *Mol Cell Endocrinol* 177:3-11.

FOR RESEARCH USE ONLY.

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