

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT S1P5 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1054-1

Lot Number: DC1054-1-031121

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human S1P5 cDNA (GenBank Accession Number: NM_030760) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Overnight seeding and thawing Media (cells will not grow and maintain function beyond): DMEM, 10% FBS

Background: The lysosphingolipid sphingosine 1-phosphate (S1P) regulates cell proliferation, apoptosis, motility, and neurite retractions. S1P and the structurally related LPA signal cells through a set of G-protein-coupled receptors known as EDG receptors. S1P5 or EDG8 is a 400-amino acid 7 transmembrane protein that shares 42-49% amino acid identity with the human S1P receptors EDG1, EDG3, and EDG5. Northern blot analysis has shown widespread expression of EDG8 in rat tissues, with strongest expression in spleen and white matter tracts of the brain.

Application: Functional assay

Figure 1

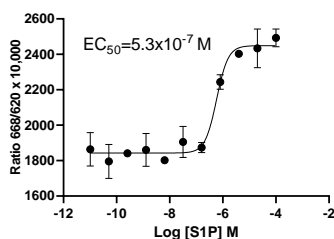


Figure 1. Dose-dependent inhibition of forskolin stimulated intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

IM *et al.* (2000) Characterization of a novel phingosine 1-phosphate receptor, Edg-8. *J Biol Chem* 275:14281-14286.

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.