

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT S1P1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DCG1047-1

Lot Number: 09/30/11

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gaqi5

Transfection: Expression vector containing full-length human S1P1 cDNA (GenBank Accession Number NM_001400) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: S1P1 (or EDG1) is a widely distributed G-protein-coupled receptor for sphingosine-1-phosphate (S1P), a blood-borne bioactive lipid. Stimulation of the S1P1 receptor triggers a G_i-linked pathway, leading to growth, survival, migration, and morphogenesis. Disruption of the S1P1 gene in mice results in embryonic lethality because of its key role within endothelial cells in regulating the coverage of blood vessels by vascular smooth muscle cells. S1P1 also mediates activation of Rac and functions as a typical chemotactic receptor.

Application: Functional assay

Figure 1

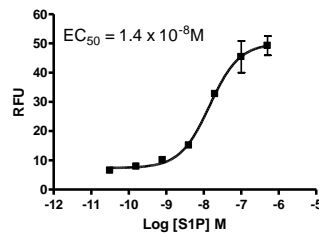


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

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.

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