

**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 ( $4 \times 10^6$ ) 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

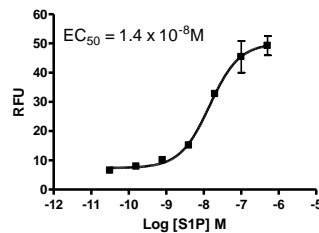
**Propagation Medium:** DMEM/F12, 10% FBS

**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<sub>i</sub>-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.