

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT S1P3 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DCG1049-1

Lot Number: CG1049-1-091610

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gaqi5

Transfection: Expression vector containing full-length human S1P3 cDNA (Genbank Accession Number NM_005226.2) 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: S1P3 (sphingosine 1-phosphate receptor 3) or EDG3 (endothelial differentiation G-protein coupled receptor 3) is a receptor for the lysosphingolipid sphingosine 1-phosphate, a bioactive lysophospholipid that elicits diverse physiological effect on most types of cells and tissues. EDG receptors mediate pro-angiogenic activities, such as endothelial cell proliferation, chemotaxis, and vessel morphogenesis. S1P3, along with S1P1, is required for directing the organization of endothelial cells into capillary networks.

Application: Functional assay

Figure 1

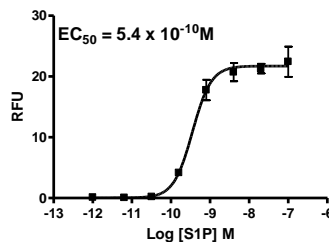


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.