

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT S1P5 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1054-1

**Lot Number:** C1054-1-082815

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**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

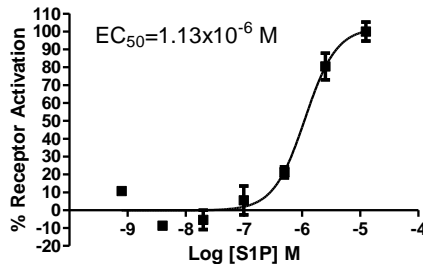
**Propagation Medium:** DMEM/F12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** Stable in culture for minimum of two months

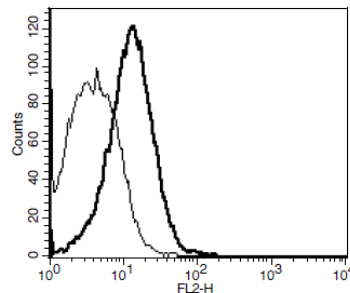
**Background:** The lysosphingolipid sphingosine 1-phosphate (S1P) regulates cell proliferation, apoptosis, motility, and neurite retractions. S1P and the structurally related LPA signal cells through a set of G-protein-coupled receptors known as EDG receptors. S1P5 or EDG8 is a 400-amino acid 7 transmembrane 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 2**



**Figure 1.** Dose-dependent inhibition of forskolin stimulated intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

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

**FOR RESEARCH USE ONLY.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.