

PRODUCT DATASHEET
ChemiScreen™ S1P₁ Lysophospholipid Membrane Preparation

CATALOG NUMBER:	HTS176M	QUANTITY:	200 units
LOT NUMBER:		VOLUME/CONCENTRATION	1 mL, 1 mg/mL

BACKGROUND: Sphingosine 1-phosphate (S1P) is a biologically active lysophospholipid that transmits signals through a family of five G-protein-coupled receptors to regulate cell proliferation, migration, cytoskeletal organization, and differentiation. S1P₁ was the first identified S1P receptor. It primarily couples to PTX-sensitive Gi/o proteins and mediates S1P-induced adenylate cyclase inhibition. Expression of S1P₁ is pervasive, including spleen, brain, heart, lung, adipose tissues, liver, thymus, kidney, and skeletal muscle (Zhang *et al.* 1999). The deletion of S1P₁ in mice results in embryonic lethality (Liu *et al.*, 2000) with death attributable to incomplete vascular maturation. Recent reports demonstrate specific roles for S1P₁ in lymphocyte recirculation/egress (Matloubian *et al.*, 2004). S1P₁ membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level of GPCR surface expression; thus, they are ideal HTS tools for screening of S1P₁ interactions with its ligands. The membrane preparation exhibits an EC₅₀ of 9.5 nM for Sphingosine 1-phosphate in a GTP_γS binding assay.

APPLICATIONS: GTP_γS Binding

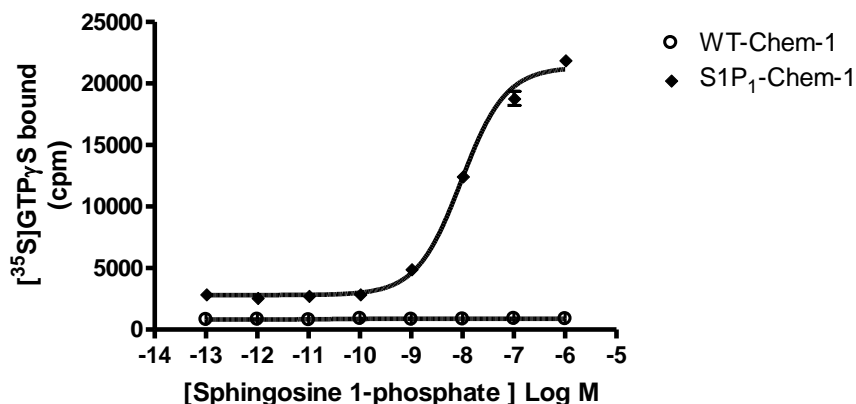


Figure 1. Binding of [³⁵S]-GTP_γS to S1P₁ membrane preparation with high GDP concentration. 5 μg/well S1P₁ Membrane Preparation (catalog # HTS176M) or Wild-Type Chem-1 Membrane Preparation (catalog # HTS000MC1) was incubated with 0.3 nM [³⁵S]-GTP_γS, 10 μM GDP and increasing amounts of unlabeled sphingosine 1-phosphate. Bound radioactivity was determined by filtration and scintillation counting.

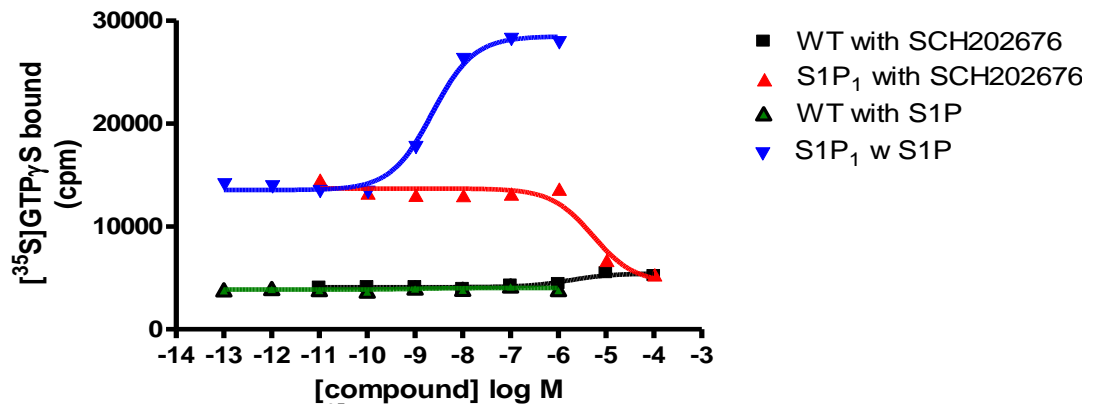


Figure 2. Binding of [³⁵S]-GTP_γS to S1P₁ membrane preparation with low GDP concentration. 5 μg/well S1P₁ Membrane Preparation (catalog # HTS176M) or Wild-Type Chem-1 Membrane Preparation (catalog # HTS000MC1) was incubated with 0.3 nM [³⁵S]-GTP_γS, 0.5 μM GDP and increasing amounts of unlabeled sphingosine 1-phosphate or SCH202676. Bound radioactivity was determined by filtration and scintillation counting. Increased constitutive activity of S1P₁ was observed with low GDP concentration, and this constitutive activity was inhibited by the broad spectrum GPCR inhibitor SCH202676.

SPECIFICATIONS: 1 unit = 5 μg

EC50 in GTP_γS binding assay by Sphingosine 1-phosphate: ~ 9.5 nM

Species: Full-length human EDG1 cDNA encoding S1P₁ (Accession Number: NM_001400)

HOST CELLS: Chem-1, an adherent cell line expressing the promiscuous G-protein, G_α15.

RECOMMENDED ASSAY CONDITIONS: Membranes are permeabilized by addition of saponin to an equal concentration by mass, then mixed with [³⁵S]-GTP_γS (final concentration of 0.3 nM) in 20 mM HEPES, pH 7.4/100 mM NaCl/10 mM MgCl₂/10 μM GDP in a nonbinding 96-well plate. Unlabeled Sphingosine 1-phosphate was added to the final concentration indicated in Figure 1 (final volume 100 μL), and incubated for 30 min at 30°C. The binding reaction is transferred to a GF/B filter plate (Millipore MAHF B1H) previously prewetted with water. The plate is washed 3 times (1 mL per well per wash) with cold 10 mM sodium phosphate, pH 7.4, then dried and counted.

Note: Performing the reaction with lower GDP concentrations (0.5 μM) results in elevated constitutive activity (Figure 2).

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 1000 cpm specific S1P-stimulated [³⁵S]-GTP_γS binding.

PRESENTATION:

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.

Packaging method: Membrane protein was adjusted to the indicated concentration in packaging buffer, rapidly frozen, and stored at -80°C.

STORAGE/HANDLING:

Store at -70°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

REFERENCES:

1. Zhang GF *et al.* (1999) Comparative analysis of three murine G-protein coupled receptors activated by sphingosine-1-phosphate. *Gene* 227: 89–99
2. Liu Y *et al.* (2000) Edg-1, the G protein–coupled receptor for sphingosine-1-phosphate, is essential for vascular maturation. *J. Clin. Invest.* 106, 951–961
3. Matloubian M *et al.* (2004) Lymphocyte egress from thymus and peripheral lymphoid organs is dependent on S1P receptor 1. *Nature* 427: 355–360

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services