



CHEMISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT S1P₂ LYSOPHOSPHOLIPID RECEPTOR

CATALOG NUMBER:	HTS078M	QUANTITY:	200 units
LOT NUMBER:	2022543	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL

BACKGROUND: Sphingosine 1-phosphate (S1P) is a bioactive lipid that binds to and activates a family of GPCRs, S1P₁₋₅ (also known as EDG receptors). Interactions between S1P and its receptors mediate cytoskeletal rearrangement and cell migration, with functional consequences in angiogenesis, lymphocyte trafficking, and smooth muscle development (Anliker and Chun, 2004). S1P₁ (Edg-1) signals exclusively through G_i, whereas S1P₂ (Edg-5) and S1P₃ (Edg-3) activate G_i, G_q and G_{12/13} (Windh *et al.*, 1999). Although S1P₁ and S1P₃ promote cell migration, S1P₂ inhibits cell migration in several cell types; these opposing functions appear to result from differences in the ability of each receptor to activate G_i (Arikawa *et al.*, 2003; Sugimoto *et al.*, 2003; Goparaju *et al.*, 2005). Studies with knockout mice indicate that S1P₂ and S1P₃ have redundant functions in maintaining vascular integrity during embryonic development (Kono *et al.*, 2004). Chemicon's S1P₂ 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 S1P₂ interactions with its ligands. The cell line exhibits EC₅₀ of 7.2nM for S1P in a GTPγS binding assay.

APPLICATIONS: GTPγS Binding and Radioligand Binding Assay.

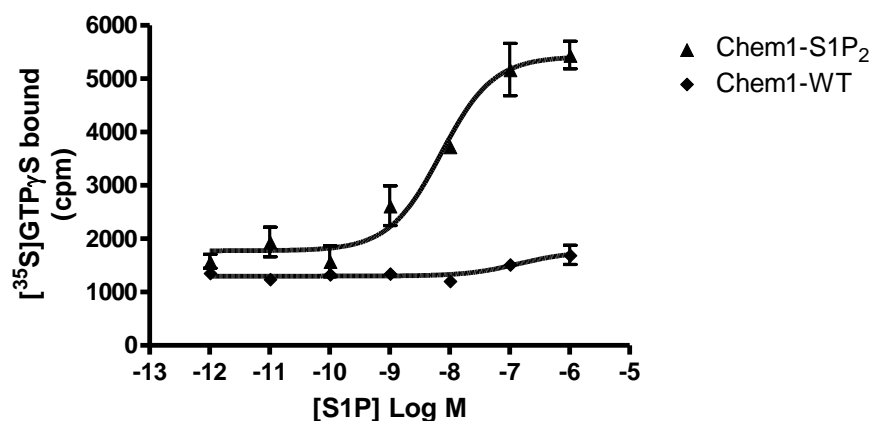


Figure 1. Binding of [³⁵S]-GTPγS to S1P₂ membrane preparation. 5 μg/well S1P₂ Membrane Preparation (catalog # HTS078M) was incubated with 0.1 nM [³⁵S]-GTPγS and increasing amounts of unlabeled S1P. Bound radioactivity was determined by filtration and scintillation counting.



SPECIFICATIONS: EC50 in GTP γ S binding assay by S1P: ~ 7.2nM

Species: human EDG-5 cDNA encoding S1P₂ (Accession Number: [NM_004230](#))

HOST CELLS: Chem-1, an adherent cell line expressing the promiscuous G-protein, G α 15.

ASSAY CONDITIONS: Membranes are permeabilized by addition of saponin to an equal concentration by mass, then mixed with [³⁵S]-GTP γ S (final concentration of 0.1 nM) in 20 mM HEPES, pH 7.4/100 mM NaCl/10 mM MgCl₂/0.5 μ M GDP in a nonbinding 96-well plate. Unlabeled S1P is added to the final concentration indicated in Figure 1 (final volume 100 μ L), 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, and washed 3 times (1 mL per well per wash) with cold 10 mM sodium phosphate, pH 7.4. The plate is dried and counted.

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

The S1P₂ membrane preparation is expected to be functional in a radioligand binding assay; however, the end user will need to determine the optimal radiolabeled ligand for use with this product.

PRESENTATION:

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol with no preservatives.
Packaging method: Membrane protein was adjusted to 1 mg/ml in packaging buffer, rapidly frozen, and stored at -80°C.

STORAGE/HANDLING:

Maintain frozen at -70°C for up to 2 years. Do not freeze and thaw.

REFERENCES:

Anliker B and Chun J (2004) Lysophospholipid G Protein-coupled Receptors. *J. Biol. Chem.* 279: 20555-20558.

Arikawa K *et al.* (2003) Ligand-dependent Inhibition of B16 Melanoma Cell Migration and Invasion via Endogenous S1P₂ G Protein-coupled Receptor. *J. Biol. Chem.* 278: 32841-32851.

Goparaju SK *et al.* (2005) The S1P₂ Receptor Negatively Regulates Platelet-Derived Growth Factor-Induced Motility and Proliferation. *Mol. Cell. Biol.* 25: 4237-4249.

Kono M *et al.* (2004) The Sphingosine-1-phosphate Receptors S1P₁, S1P₂, and S1P₃ Function Coordinately during Embryonic Angiogenesis. *J. Biol. Chem.* 279: 29367-29373

Sugimoto N *et al.* (2003) Inhibitory and Stimulatory Regulation of Rac and Cell Motility by the G_{12/13}-Rho and G_i Pathways Integrated Downstream of a Single G Protein-Coupled Sphingosine-1-Phosphate Receptor Isoform. *Mol. Cell. Biol.* 23: 1534-1545.

Windh RT *et al.* (1999) Differential Coupling of the Sphingosine 1-Phosphate Receptors Edg-1, Edg-3, and H218/Edg-5 to the G_i, G_q, and G₁₂ Families of Heterotrimeric G Proteins. *J. Biol. Chem.* 274: 27351-27358.



Important Note: *During shipment, small volumes of product will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ L or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.*

**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.

©2002 - 2013: Merck KGaA, Darmstadt. All rights reserved. No part of these works may be reproduced in any form without permission in writing.