

Product Information

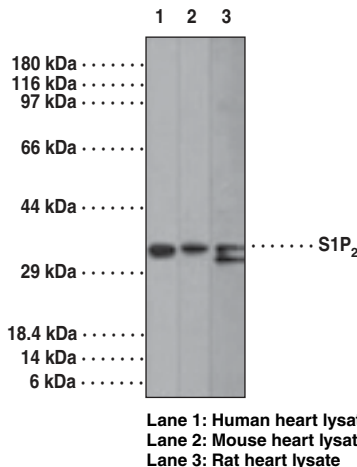


S1P₂ Polyclonal Antibody

Item No. 13488

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Synonyms:** EDG-5, S1PR2, Sphingosine-1-phosphate Receptor 2
- Antigen:** Synthetic peptide from human S1P₂ amino acids 325-343 (LLPLRSSSLERGMHMPTS)
- Host:** Rabbit
- Cross Reactivity:** (+) Human, monkey, mouse, and rat S1P₂
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 1-2 µg/ml.

S1P₂ is a membrane protein belonging to the EDG family of extracellular G protein-coupled receptors. It acts as a receptor for the bioactive lipid sphingosine 1-phosphate (S1P) and participates in S1P-induced cell proliferation, survival, and transcriptional activation. It shares 92% amino acid identity with rat H218 and approximately 44% identity with the human S1P receptors EDG-1 and EDG-3. Though the role of S1P₂ in vascular permeability is not clear, reports suggest that it regulates PTEN by a Rho GTPase-dependent pathway to inhibit cell migration. S1P₂ also plays an important role in normal vascularization and hypoxia-triggered pathological angiogenesis of the mouse retina. The gene for human S1P₂ protein is localized on chromosomal region 19p13.2. The protein is usually expressed in human brain and heart.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13488

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/14/2011