

Product Information

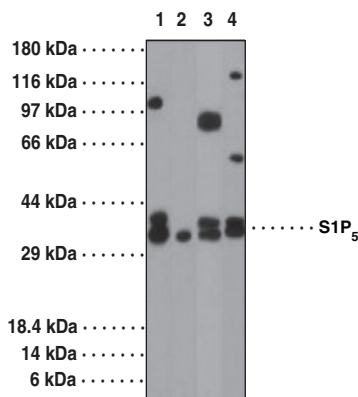


S1P₅ Polyclonal Antibody

Catalog No. 13490

- 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-8, S1PR5, Sphingosine-1-phosphate Receptor 5
- Antigen:** Synthetic peptide from human S1P₅ within the region of amino acids 1-50
- Host:** Rabbit
- Cross Reactivity:** (+) Canine, human, monkey, murine, and rat S1P₅
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 1-3 µg/ml and 5-10 µg/ml in immunohistochemistry (paraffin).

Sphingosine 1-phosphate (S1P) receptor EDG-8 is a multi-pass membrane protein belonging to the G protein-coupled receptor 1 family and coupled to both the G_{i/oα} and G₁₂ subclass of heteromeric G-proteins. EDG-8 acts as a receptor for the bioactive sphingolipid S1P as well as for lysophosphatidic acid (LPA) and gives diverse physiological effects on most types of cells and tissues. EDG-8 may also play a regulatory role in the transformation of radial glial cells into astrocytes and affect proliferative activity of these cells. It is widely expressed in white matter tracts of brain, spleen, peripheral blood leukocytes, placenta, lung, aorta, and fetal spleen. Over expression of EDG-8 leads to a lymphoproliferative disorder often associated with the autoimmune disease known as leukemic large granular lymphocyte (LGL).



Lane 1: Human brain lysate (Absence of immunizing peptide)
Lane 2: Human brain lysate (Presence of immunizing peptide)
Lane 3: Murine S1P₅
Lane 4: Rat S1P₅

Related Products

S1P₂ Polyclonal Antibody - Cat. No. 13488 • S1P₄ Peptide Affinity-Purified Polyclonal Antibody - Cat. No. 13489 • LPA₃ Polyclonal Antibody - Cat. No. 10004840 • S1P₁ Polyclonal Antibody - Cat. No. 10005228 • LPA₁ Polyclonal Antibody - Cat. No. 10005280 • S1P₃ Polyclonal Antibody - Cat. No. 10006373

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, 10/19/2009