

Product Information

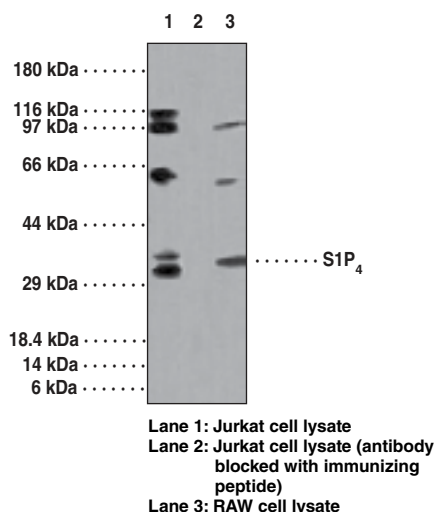


S1P₄ Polyclonal Antibody

Item No. 13489

Contents:	This vial contains 100 µg of peptide affinity-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonyms:	EDG-6, S1PR4, Sphingosine-1-phosphate Receptor 4
Antigen:	Synthetic peptide from human S1P ₄ within the region of amino acids 1-50
Host:	Rabbit
Cross Reactivity:	(+) Human and mouse S1P ₄
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	The recommended starting dilution for western blot is 3.0-5.0 µg/ml and 5 µg/ml in IF and immunocytochemistry.

EDG-6 is a member of the endothelial differentiation gene (EDG) G protein-coupled receptor family and is a receptor for sphingosine 1-phosphate (S1P). Binding of S1P to EDG-6 activates Erk, leading to the activation of the transcription factor Elk1 and thus the mitogen-activated protein kinase (MAPK) signal transduction pathway. The protein is specifically expressed in fetal and adult lymphoid and hematopoietic tissue as well as in lung. Human EDG-6 gene is intronless and is mapped to chromosome 19p13.3 at the D19S120 marker.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13489

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, 12/15/2011