

Product Information



Serine Palmitoyltransferase Blocking Peptide

Item No. 10006620

Sphingolipids play essential roles in various cellular events, including proliferation, differentiation, senescence, apoptosis, and inflammatory responses.¹ Serine palmitoyltransferase (SPT) is the initial and rate-limiting enzyme in the *de novo* sphingolipid biosynthesis, and thus regulates the level of sphingolipids in cells.² An immunohistochemical study revealed widespread distribution of the enzyme with pronounced expression in brain and digestive tract.³ Two subunits, SPT1 and SPT2 at a stoichiometry of 1:1, are involved in the enzymatic activity of SPT.⁴ Cayman Chemical's serine palmitoyltransferase polyclonal antibody recognizes SPT2, the long chain subunit of the enzyme. The antibody stains mainly cell nuclei and both cell nuclei and cytoplasm of a subpopulation of cells in formalin-fixed, paraffin-embedded rat brain tissue. The nuclear localization of SPT2 may suggest that SPT2 associates with another nuclear protein or is modified and transported to the nucleus.²

Laboratory Procedures

The SPT Blocking Peptide (human serine palmitoyltransferase subunit SPT2 amino acids 548-562) can be used in conjunction with Cayman's SPT Polyclonal Antibody (Item No. 10005260) to block protein-antibody complex formation during immunochemical analysis of SPT.

Store this peptide solution at -20°C. It will be stable for at least two years. To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the SPT Polyclonal Antibody (Item No. 10005260) and blocking peptide together in a 1:2 (v/v) ratio in a microfuge tube. For example, mix 40 µl of antibody and 80 µl of peptide.*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

References

1. Mathias, S., Peña, L.A., and Kolesnick, R.N. Signal transduction of stress *via* ceramide. *Biochem J.* **335**, 465-480 (1998).
2. Carton, J.M., Uhlinger, D.J., Batheja, A.D., *et al.* Enhanced serine palmitoyltransferase expression in proliferating fibroblasts, transformed cell lines, and human tumors. *J. Histochem. Cytochem.* **51(6)**, 715-726 (2003).
3. Batheja, A.D., Uhlinger, D.J., Carton, J.M., *et al.* Characterization of serine palmitoyltransferase in normal human tissues. *J. Histochem. Cytochem.* **51(5)**, 687-696 (2003).
4. Hanada, K., Hara, T., and Nishijima, M. Purification of the serine palmitoyltransferase complex responsible for sphingoid base synthesis by using affinity peptide chromatography techniques. *J. Biol. Chem.* **275(12)**, 8409-8415 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10006620

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, 07/19/2012

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