

Product Information



ACAT-1 Blocking Peptide

Catalog No. 10005090

Acyl-coenzyme A: cholesterol acyltransferase-1 (ACAT-1) catalyzes the formation of cholesterol esters from cholesterol and long chain fatty acyl-coenzyme A. ACAT-1 may be involved in maintaining appropriate membrane free cholesterol levels and in lipid droplet formation, and thus play a role in the development of atherosclerosis.^{1,2} ACAT-1 is ubiquitously expressed, with highest levels observed in sebaceous glands, steroidogenic tissues, and macrophages.³⁻⁵ This intracellular enzyme is located at the endoplasmic reticulum and has as many as eight transmembrane domains.¹ Human ACAT-1 has 550 amino acids with an estimated molecular weight of 65 kDa.⁵ Cayman's affinity purified antibody recognizes a 50 kDa band in most cultured cell lines as reported in the literature.^{4,6,7} In addition, the antibody reacts with a native form of ACAT-1 at about 65 kDa in liver tissues.

Laboratory Procedures

This vial contains 200 µg peptide in 200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. Store aliquots at -20°C; they will be good for up to two years. The ACAT-1 blocking peptide (human ACAT-1 amino acids 6-23) can be used in conjunction with Cayman's ACAT-1 Polyclonal Antibody (Catalog No. 100028) to block protein-antibody complex formation during immunochemical analysis of ACAT-1

Reconstitute the lyophilized peptide with 200 µl of PBS or distilled water. 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 ACAT-1 Polyclonal Antibody (Catalog No. 100028) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 20 µl of peptide.*
2. Incubate for 1 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. Rudel, L.L., Lee, R.G., and Cockman, T.L. Acyl coenzyme A: cholesterol acyltransferase types 1 and 2: structure and function in atherosclerosis. *Curr. Opin. Lipidol.* **12**, 121-127 (2001).
2. Linton, M.F. and Fazio, S. Macrophages, inflammation, and atherosclerosis. *Int. J. Obes.* **27**, S35-S40 (2003).
3. Namatame, I., Tomoda, H., Ishibashi, S., *et al.* Antiatherogenic activity of fungal beauveriolides, inhibitors of lipid droplet accumulation in macrophages. *Proc. Natl. Acad. Sci. USA* **101**(3), 737-742 (2004).
4. Lee, R.G., Willingham, M.C., Davis, M.A., *et al.* Differential expression of ACAT1 and ACAT2 among cells within liver, intestine, kidney, and adrenal of nonhuman primates. *J. Lipid Res.* **41**, 1991-2001 (2000).
5. Chang, C.C.Y., Huh, H.Y., Cadigan, K.M., *et al.* Molecular cloning and functional expression of human acyl-coenzyme A: Cholesterol acyltransferase cDNA in mutant chinese hamster ovary cells. *J. Biol. Chem.* **268**(28), 20747-20755 (1993).
6. Chang, C.C.Y., Lee, C.-Y.G., Chang, E.T., *et al.* Recombinant acyl-CoA: cholesterol acyltransferase-1 (ACAT-1) purified to essential homogeneity utilizes cholesterol in mixed micelles or in vesicles in a highly cooperative manner. *J. Biol. Chem.* **273**(52), 35132-35141 (1998).
7. Chang, C.C.Y., Chen, J., Thomas, M.A., *et al.* Regulation and immunolocalization of acyl-coenzyme A: cholesterol acyltransferase in mammalian cells as studies with specific antibodies. *J. Biol. Chem.* **270**(49), 29532-29540 (1995).

Related Product

ACAT-1 Polyclonal Antibody - Cat. No. 100028

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSDS supervisor at 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, 11/03/2007

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