

Product Information



sPLA₂ (human Type IIA) Affinity Sorbent

Item No. 485009

IMPORTANT

Water used to prepare all EIA reagents and buffers must be deionized and free of trace organic contaminants ('UltraPure'). Use activated carbon filter cartridges or other organic scavengers. Glass distilled water (even if double distilled), HPLC-grade water, and sterile water (for injections) are not adequate for EIA. UltraPure water may also be purchased (Item No. 400000).

Instructions for Use

This vial contains 1 ml of an sPLA₂ (human synovial) affinity sorbent (Mouse Anti-sPLA₂ covalently bound to Sepharose 4B). One milliliter of sorbent is capable of binding approximately 4 µg sPLA₂. The antibody was raised against type IIA sPLA₂, but may also bind other sPLA₂s.

The sorbent may be used for the removal of sPLA₂ from various media. Because Sepharose 4B is a gel filtration matrix, it is not advisable to use the sorbent in a column format to remove sPLA₂ from small volumes of sample since other small molecules may also be removed. Instead, add the sorbent to the sample and vortex. Centrifuge at 500 x g to precipitate the sorbent and pour off the supernatant, which should now be free of sPLA₂. To regenerate the sorbent simply wash with 10 ml UltraPure water, 10 ml 50 mM glycine • HCl, pH 2.5, containing 0.1% Triton X-100 and 0.15 M NaCl, or 50 mM Triethanolamine, pH 11.5, containing 0.1% Triton X-100 and 0.15 M NaCl, and 10 ml water. Using a small fitted filter will speed this process.

Related Products

sPLA₂ (human Type IIA) Monoclonal Antibody - Item No. 160500 • sPLA₂ (human Type IIA) Polyclonal Antiserum - Item No. 160502 • sPLA₂ (human Type IIA) EIA Kit - Item No. 585000 • sPLA₂ Assay Kit - Item No. 765001

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, 07/29/2010

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