

# Product Information



## SPE Cartridges (Mixed Bed)

Item No. 400021

The SPE Cartridges (mixed bed) are mixed-mode cartridges packed with non-polar and strong anion exchange chemistries. As this product relates to Cayman's product line, this column has been utilized for the purification of 11-dehydro thromboxane B<sub>2</sub> (11-dehydro TXB<sub>2</sub>) from urine samples.<sup>1</sup> If your samples are being purified and analyzed for compounds other than 11-dehydro TXB<sub>2</sub>, you may want to contact Cayman directly for appropriate technical information.

### Purification Protocol

Solutions needed:

1. Buffer A: 63 mM ammonium bicarbonate, pH 8.6
2. Buffer B: 1:100 dilution of Buffer A (do not adjust pH of diluted buffer)
3. Methanol
4. Elution Solution: methanol containing 2% formic acid; made fresh just before use
5. 11-dehydro Thromboxane B<sub>2</sub> EIA Buffer Concentrate (10X) (Item No. 419507); dilute as instructed on the insert

Procedure (based on 1 ml of sample):

1. Dilute centrifuged samples with an equal volume of Buffer A and incubate at room temperature for three hours.
2. Condition column with 3 ml methanol, followed by 3 ml deionized water. Do not allow the column to go dry.
3. Apply the sample to the column using gravity flow.
4. Rinse the column sequentially with 3 ml of Buffer B, 3 ml deionized water, and then with 6 ml methanol.
5. Elute the 11-dehydro TXB<sub>2</sub> using 2 ml of elution solution *NOTE: It is critical that the elution solution be used within 1-2 hours of preparation.*
6. Evaporate the sample to dryness under a gentle stream of dry nitrogen.
7. Resuspend the sample in 1 ml of diluted 11-dehydro TXB<sub>2</sub> EIA Buffer.

### Reference

1. Perneby, C., Granstrom, E., Beck, O., *et al.* Optimization of an enzyme immunoassay for 11-dehydro-thromboxane B<sub>2</sub> in urine: Comparison with GC-MS. *Thromb. Res.* **96**, 427-436 (1999).

### Related Products

SPE Cartridges (C-18) - Item No. 400020 • 11-dehydro Thromboxane B<sub>2</sub> EIA Kit - Item No. 519501

**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, 10/18/2010

## 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