

Product Information



Prostaglandin E₂ Affinity Purification Kit (4 ml)

Item No. 500450

Table of Contents

Item No.	Item Name	1 ea Quantity/Size	5 ea Quantity/Size
400220	Eicosanoid Affinity Column Buffer (5X)	1 vial/50 ml	1 vial/250 ml
400230	Eicosanoid Affinity Column Elution Solution	1 vial/50 ml	1 vial/250 ml
400056	Prostaglandin E ₂ Affinity Column (4 ml)	1 column/0.5 ml sorbent in a 4 ml column	5 columns/0.5 ml sorbent in a 4 ml column

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 975-3999. We cannot accept any returns without prior authorization.

Precautions

Please read these instructions carefully before beginning the purification.

For research use only. Not for human or diagnostic use.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888

Fax: 734-971-3641

Email: techserv@caymanchem.com

Hours: M-F 8:00 AM to 5:30 PM EST

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored at 4°C and used before the expiration date indicated on the outside of the box.

Reagent Preparation

1. Eicosanoid Affinity Column Buffer (5X) (Item No. 400220)

Dilute 1:5 with UltraPure water (*i.e.*, dilute 50 ml vials to a total volume of 250 ml and the 250 ml vials to a total volume of 1,250 ml).

2. Prostaglandin E₂ Affinity Column (Item No. 400056)

Remove the top stopper and then gently remove the bottom plug. Be certain to remove the top stopper first to avoid air bubbles being drawn into the packing material. Allow the storage solution to pass through the gel and then wash two times with 2 ml of Column Buffer. The column is now ready to use.

3. Eicosanoid Affinity Column Elution Solution (Item No. 400230)

This solution is ready to use.

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, 10/26/2011

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

Product Information



4. EIA Buffer

Prepare a 1.0 M phosphate buffer solution by combining 133 g K_2HPO_4 and 32.15 g KH_2PO_4 and diluting to a total volume of 1.0 liter with UltraPure water. The pH of this solution will be 7.4. To 100 ml of this buffer, add the following: 100 mg sodium azide, 23.4 g sodium chloride, 370 mg tetrasodium EDTA, and 1 g bovine serum albumin (Sigma A7030 or equivalent). Stir at room temperature until completely dissolved and dilute to a total volume of 1.0 liter with UltraPure water; the pH of this solution will be ~7.0. This buffer may also be purchased as a 10X concentrated buffer (Item No. 400060).

Purification Protocol

1. All samples must be free of particulates and precipitates. This can be achieved by either filtration or centrifugation. All samples must be at approximately neutral pH (6.5-7.5). Urine samples should be centrifuged briefly to remove sediment and may be applied directly to the column. Plasma samples should be diluted 1:5 with Column Buffer and applied to the column. Allow the entire sample to pass through the packing material. The binding capacity of the column is 5 ng of prostaglandin E_2 (PGE_2). Your sample should contain an amount of PGE_2 equal to 20-50% of the column capacity.
2. Wash the column with 2 ml of Column Buffer, followed by 2 ml UltraPure water. Allow all of the water to pass through the packing material. Discard both of these washes.
3. Place a test tube capable of holding at least 5 ml under the column. Elute the PGE_2 from the column by adding 5 ml of Elution Solution and allowing it to pass through the packing material.
4. Evaporate the Elution Solution to dryness either by vacuum centrifugation or by evaporation under a stream of dry nitrogen. It is imperative that all of the organic solvent be removed as even trace quantities may adversely affect any immunoassay.
5. Immediately dissolve the samples in the buffer or solvent appropriate for your application. If you are assaying the sample with any of our PGE_2 EIA Kits (Item No. 514010 or 500141), dissolve the sample in EIA Buffer. The amount of EIA Buffer to use depends on the original sample volume and the expected concentration of PGE_2 in the sample. A dilute sample may be concentrated by dissolving the residue in a smaller volume of EIA Buffer than the original sample volume.

Recovery

Urine and plasma recoveries average >90% with a variance of <20%. If your application requires greater precision, samples may be spiked with tritiated PGE_2 prior to purification. For a protocol on recovery estimation using tritiated PGE_2 , please see the PGE_2 EIA Kit-Monoclonal booklet (Item No. 514010).

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

Product Information



Specificity

Prostaglandin E ₂	100%	tetranor-PGEM	<0.01%
Prostaglandin E ₂ Ethanolamide	100%	tetranor-PGFM	<0.01%
Prostaglandin E ₂ -1-glyceryl ester	100%	Prostaglandin A ₁	<0.01%
Prostaglandin E ₃	43%	8- <i>iso</i> Prostaglandin A ₂	<0.01%
Prostaglandin E ₁	18.7%	Prostaglandin A ₃	<0.01%
8- <i>iso</i> Prostaglandin E ₂	2.5%	Prostaglandin B ₁	<0.01%
Sulprostone	1.25%	Prostaglandin B ₂	<0.01%
6-keto Prostaglandin F _{1α}	1%	Prostaglandin D ₂	<0.01%
8- <i>iso</i> Prostaglandin F _{2α}	0.25%	11-deoxy Prostaglandin E ₂	<0.01%
Prostaglandin A ₂	0.04%	16,16-dimethyl Prostaglandin E ₂	<0.01%
13,14-dihydro-15-keto Prostaglandin E ₂	0.02%	19(R)-hydroxy Prostaglandin E ₂	<0.01%
Arachidonic acid	<0.01%	20-hydroxy Prostaglandin E ₂	<0.01%
Arachidonoyl Ethanolamide	<0.01%	15-keto Prostaglandin E ₂	<0.01%
O-Arachidonoyl Ethanolamide	<0.01%	Prostaglandin F _{1α}	<0.01%
Butaprost	<0.01%	Prostaglandin F _{2α}	<0.01%
10(E),12(Z)-Conjugated Linoleic Acid	<0.01%	13,14-dihydro-15-keto Prostaglandin F _{2α}	<0.01%
8(S),15(S)-DiHETE	<0.01%	Prostaglandin F _{3α}	<0.01%
Leukotriene B ₄	<0.01%	Prostaglandin J ₂	<0.01%
Misoprostol	<0.01%	15-deoxy-Δ ^{12,14} -Prostaglandin J ₂	<0.01%
Misoprostol (free acid)	<0.01%	Thromboxane B ₂	<0.01%

Related Products

Eicosanoid Affinity Column Buffer (5X) - Item No. 400220
Eicosanoid Affinity Column Elution Solution - Item No. 400230
Prostaglandin E₂ - Item No. 14010
Prostaglandin E₂ Affinity Column (4 ml) - Item No. 400056
Prostaglandin E₂ Affinity Column (20 ml) - Item No. 400057
Prostaglandin E₂ Affinity Purification Kit (20 ml) - Item No. 514018
Prostaglandin E₂ Affinity Sorbent - Item No. 414020
Prostaglandin E₂ EIA Kit - Item No. 514010
Prostaglandin E₂ Express EIA Kit - Item No. 500141
UltraPure Water - Item No. 400000

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