

Product Information



Prostaglandin Metabolite HPLC Mixture

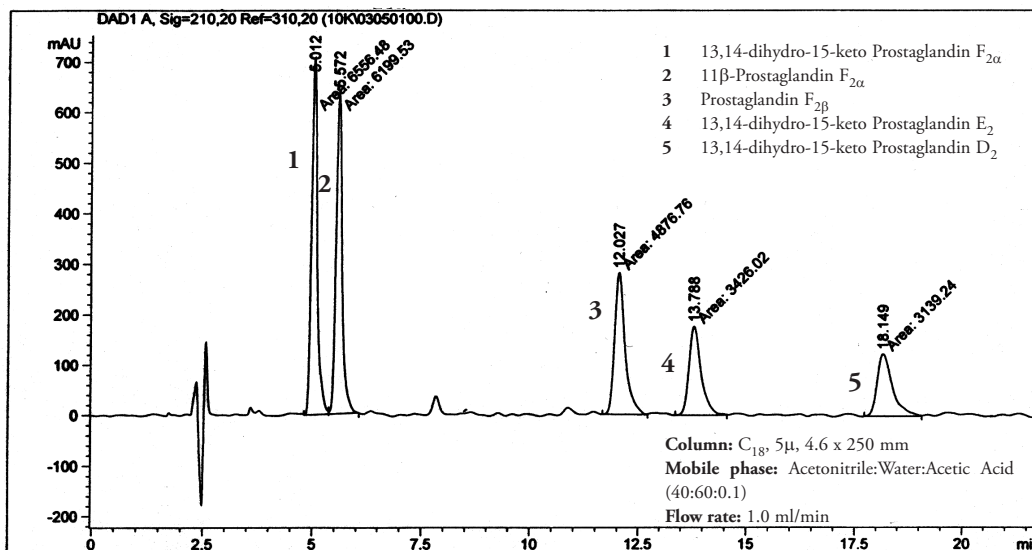
Catalog No. 10005

Purity: ≥98% for each compound
Stability: ≥6 months at -20°C
Supplied as: A solution in methyl acetate (100 µg of each compound)

Laboratory Procedures

The Prostaglandin Metabolite HPLC Mixture (PG metabolite HPLC mixture) is supplied as a solution in methyl acetate. We suggest that they be stored as supplied at -20°C. It should be stable for at least six months.

This HPLC mixture contains the primary metabolites of five commonly encountered cyclooxygenase products. In primates, the lung is an active site of PG inactivation *via* 15-hydroxy PG dehydrogenase (15-hydroxy PGDH).¹ E and F-series PGs are good substrates for 15-hydroxy PGDH, which oxidizes the C-15 hydroxyl group to give the corresponding 15-keto compounds.² This oxidation is followed rapidly by Δ^{13} -reductase removal of the 13,14-double bond, to give 13,14-dihydro-15-keto PG metabolites.³ While the primary PGs are often undetectable in plasma due to their rapid inactivation, the 13,14-dihydro-15-keto metabolites can be present at concentrations of several hundred pg/ml.⁴ In addition, 9- and 11-keto PG-reductases have been reported, especially in the case of PGD₂, which produce F-series PG metabolites with a β -hydroxyl group on the ring.⁵ These metabolites span a wide polarity range and separate well on reversed phase HPLC.



References

1. Hansen, H.S. *Prostaglandins* 12, 647-679 (1976).
2. Sun, F.F., Armour, S.B., Bockstanz, V.R., *et al. Adv. Prostaglandin Thromboxane Res.* 1, 163-169 (1976).
3. Basu, S. and Stjernschantz, J. *Prostaglandins Leukot. Essent. Fatty Acids* 57, 305-310 (1997).
4. Granström, E., Hamberg, M., Hansson, G., *et al. Prostaglandins* 19, 933-945 (1980).
5. Ellis, C.K., Smigel, M.D., Oates, J.A., *et al. J. Biol. Chem.* 254, 4152-4163 (1979).

Related Products

Cyclopentenone Prostaglandin HPLC Mixture - Cat. No. 10000 • Primary Eicosanoid HPLC Mixture - Cat. No. 10001 • Prostaglandin HPLC Mixture - Cat. No. 10002 • Vasoactive Eicosanoid HPLC Mixture - Cat. No. 10003 • 13,14-dihydro-15-keto Prostaglandin D₂ - Cat. No. 12610 • 13,14-dihydro-15-keto Prostaglandin E₂ - Cat. No. 14650 • Prostaglandin F_{2β} - Cat. No. 16410 • 11β-Prostaglandin F_{2α} - Cat. No. 16520 • 13,14-dihydro-15-keto Prostaglandin F_{2α} - Cat. No. 16670

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 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, 08/04/2009

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