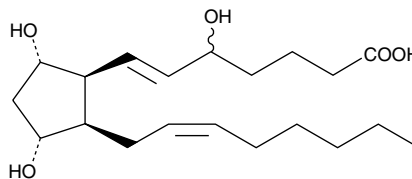


Product Information



5-*i*PF_{2α}-VI
Item No. 16300

CAS Registry No.: 180469-63-0
Formal Name: 5,9α,11α-trihydroxy-(8β)-prosta-6E,14Z-dien-1-oic acid
Synonym: *i*PF_{2α}-I
MF: C₂₀H₃₄O₅
FW: 354.5
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that (5R,S)-*i*PF_{2α}-VI be stored as supplied at -20°C. It should be stable for at least one year.

(5R,S)-*i*PF_{2α}-VI is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of (5R,S)-*i*PF_{2α}-VI in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of (5R,S)-*i*PF_{2α}-VI is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (5R,S)-*i*PF_{2α}-VI in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Isoprostanes are prostaglandin (PG)-like products of free-radical induced lipid peroxidation.¹ Although the isoprostanes derived from arachidonic acid are the best characterized, many other polyunsaturated fatty acids can form isoprostanes.² (5R,S)-*i*PF_{2α}-VI is one of dozens of possible stereo- and regioisomeric isoprostanes which can be formed from arachidonic acid. To date, the most extensively studied of these is 8-isoprostane (8-*epi*-PGF_{2α}, *i*PF_{2α}-III).^{3,4} However, 8-isoprostane is a minor isoprostane constituent when compared to some of the other isomers which form in natural conditions of oxidative stress.⁵ (5R,S)-*i*PF_{2α}-VI is an isoprostane from the unique Type VI class of isoprostanes. This class has been shown to be one of the major isoprostane products, in contrast to 8-isoprostane. In addition to being produced in greater abundance than 8-isoprostane, Type VI isoprostanes form internal lactones, which facilitates their extraction and purification from biological samples.⁵⁻⁸

References

1. Morrow, J.D., Hill, K.E., Burk, R.F., *et al. Proc. Natl. Acad. Sci. USA* **87**, 9383-9387 (1990).
2. Parchmann, S. and Mueller, M.J. *J. Biol. Chem.* **273**, 32650-32655 (1998).
3. Delanty, N., Reilly, M., Pratico, D., *et al. Br. J. Clin. Pharmacol.* **42**, 15-19 (1996).
4. Reilly, M.P., Barry, P., Lawson, J.A., *et al. Fibrinolysis & Proteolysis* **11**, 81-84 (1997).
5. Reilly, M.P., Pratico, D., Delanty, N., *et al. Circulation* **98**, 2822-2828 (1998).
6. Li, H., Lawson, J.A., Reilly, M., *et al. Proc. Natl. Acad. Sci. USA* **96(23)**, 13381-13386 (1999).
7. Lawson, J.A., Li, H., Rokach, J., *et al. J. Biol. Chem.* **273**, 29295-29301 (1998).
8. Praticò, D., Barry, O.P., Lawson, J.A., *et al. Proc. Natl. Acad. Sci. USA* **95**, 3449-3454 (1998).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16300

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, 11/04/2011

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