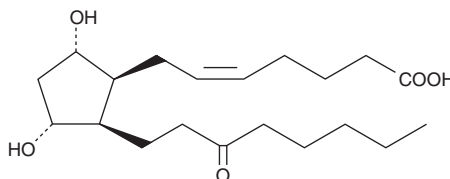


Product Information



8-*iso*-13,14-dihydro-15-keto Prostaglandin F_{2α} Catalog No. 16380

Cas Registry No.: 191919-02-5
Formal Name: 9α,11α-dihydroxy-15-oxo-(8β)-prost-5Z-en-1-oic acid
MF: C₂₀H₃₄O₅
FW: 354.5
Purity: ≥98%
Stability: ≥1 year at -80°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 8-*iso*-13,14-dihydro-15-keto Prostaglandin F_{2α} (8-*iso*-13,14-dihydro-15-keto PGF_{2α}) be stored as supplied at -80°C. It will be stable for at least one year.

8-*iso*-13,14-dihydro-15-keto PGF_{2α} is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, or dimethyl formamide purged with an inert gas can be used. The solubility of 8-*iso*-13,14-dihydro-15-keto PGF_{2α} in these solvents is approximately 100 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 aqueous solution of 8-*iso*-13,14-dihydro-15-keto PGF_{2α} is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 8-*iso*-13,14-dihydro-15-keto PGF_{2α} in PBS (pH 7.2) is approximately 10 mg/ml. For greater aqueous solubility, 8-*iso*-13,14-dihydro-15-keto PGF_{2α} can be directly dissolved in 10 mM Na₂CO₃ (6.5 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. Store aqueous solutions of 8-*iso*-13,14-dihydro-15-keto PGF_{2α} on ice and use within 12 hours of preparation. Although the aqueous solutions of 8-*iso*-13,14-dihydro-15-keto PGF_{2α} may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

8-*iso*-13,14-dihydro-15-keto PGF_{2α} is a metabolite of the isoprostane, 8-*iso* PGF_{2α} (8-isoprostane) in rabbits, monkeys, and humans.¹ 8-*iso* PGF_{2α} is a PG-like product of non-specific lipid peroxidation.² In both humans and monkeys, exogenously infused 8-isoprostane is converted primarily to metabolites having 2 or 4 carbon atoms removed from the top side chain by β-oxidation.¹ A similar pattern is observed when tritiated 8-isoprostane is infused into rabbits.³ Early in the infusion (within 10 minutes) 8-*iso*-13,14-dihydro-15-keto PGF_{2α} was a significant component of the metabolite profile, which was comprised mostly of dinor 8-isoprostane metabolites. 8-*iso*-13,14-dihydro-15-keto PGF_{2α} weakly inhibits the U-46619 or collagen-induced aggregation of human platelets. A number of the E-series isoprostanes are much more potent inhibitors of platelet aggregation than 8-*iso*-13,14-dihydro-15-keto PGF_{2α}.⁴

References

1. Chiabrando, C., Valagussa, A., Rivalta, C., *et al.* Identification and measurement of endogenous β-oxidation metabolites of 8-*epi*-prostaglandin F_{2α}. *J. Biol. Chem.* **274**, 1313-1319 (1999).
2. Morrow, J.D., Hill, K.E., Burk, R.F., *et al.* A series of prostaglandin F₂-like compounds are produced *in vivo* in humans by a non-cyclooxygenase, free radical-catalyzed mechanism. *Proc. Natl. Acad. Sci. USA* **87**, 9383-9387 (1990).
3. Basu, S. Metabolism of 8-*iso*-prostaglandin F_{2α}. *FEBS Lett.* **428**, 32-36 (1998).
4. Cranshaw, J.H., Evans, T.W., and Mitchell, J.A. Characterization of the effects of isoprostanes on platelet aggregation in human whole blood. *Br. J. Pharmacol.* **132**, 1699-1706 (2001).

Related Products

Prostaglandin F_{2α} - Cat. No. 16010 • 8-*iso* Prostaglandin F_{2α} - Cat. No. 16350 • 8-*iso*-13,14-dihydro-15-keto Prostaglandin F_{2α} - Cat. No. 16380 • 8-*iso*-15-keto Prostaglandin F_{2α} - Cat. No. 16390 • 13,14-dihydro-15-keto Prostaglandin F_{2α} - Cat. No. 16670 • 15-hydroxy Prostaglandin Dehydrogenase Polyclonal Antibody - Cat. No. 160615

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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 MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

Copyright Cayman Chemical Company, 03/19/2008

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