

Product Information



8,12-*iso*-iPF_{2α}-VI-1,5-lactone

Item No. 10312

Formal Name: 6-((E)-2-((1R,2S,3R,5S)-3,5-dihydroxy-2-((Z)-oct-2-enyl)cyclopentyl)vinyl)tetrahydro-2H-pyran-2-one

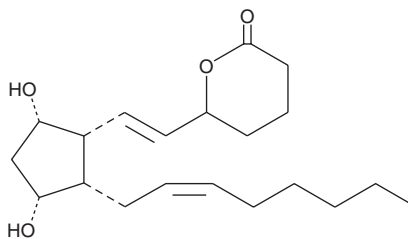
MF: C₂₀H₃₂O₄

FW: 336.2

Purity: ≥98%

Stability: ≥1 year at -80°C

Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 8,12-*iso*-iPF_{2α}-VI-1,5- lactone be stored as supplied at -80°C. It should be stable for at least one year.

8,12-*iso*-iPF_{2α}-VI-1,5- lactone 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, and dimethyl formamide purged with an inert gas can be used. The solubility of 8,12-*iso*-iPF_{2α}-VI-1,5- lactone 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 solution of 8,12-*iso*-iPF_{2α}-VI-1,5- lactone is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 8,12-*iso*-iPF_{2α}-VI-1,5- lactone in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

F₂ isoprostanes (F₂-iPs) are thought to arise from the free radical-mediated peroxidation of phospholipid-bound arachidonic acid.¹ They are cleaved, presumably by an unidentified phospholipase A₂, and are found in the circulation and the urine. 8,12-*iso*-iPF_{2α}-VI-1,5-lactone is a racemic mixture of the lactone form of the free acid, 8,12-*iso*-iPF_{2α}-VI. Previously called iPF_{2α}-I, the free acid form, iPF_{2α}-VI, is the most abundant F₂-iP regioisomer measured in the urine of rats treated with CCl₄ to induce lipid peroxidation.^{2,3} iPF_{2α}-VI is the only regioisomer that undergoes lactonization, and this occurs slowly *in vivo* or can be driven chemically.⁴ The less polar lactone is readily separated from the free acid forms of iPF_{2α}.⁴ While the level of iPF_{2α}-VI in plasma, urine, and organs is used as a biomarker for oxidative stress, some F₂-iPs also evoke significant biological effects.⁵⁻⁷ It is not known if 8,12-*iso*-iPF_{2α}-VI-1,5-lactone has important physiological effects.

References

1. Lawson, J.A., Li, H., Rokach, J., *et al.* *J. Biol. Chem.* **273**, 29295-29301 (1998).
2. Morrow, J.D., Hill, K.E., Burk, R.F., *et al.* *Proc. Natl. Acad. Sci. USA* **87**, 9383-9387 (1990).
3. Waugh, R.J., Morrow, J.D., Roberts, L.J., II, *et al.* *Free Radic. Biol. Med.* **23**, 943-954 (1997).
4. Adiyaman, M., Lawson, J.A., Khanapure, S.P., *et al.* *Anal. Biochem.* **262**, 45-56 (1998).
5. Practico, D., Barry, O.P., Lawson, J.A., *et al.* *Proc. Natl. Acad. Sci. USA* **95**, 3449-3454 (1998).
6. Van Eck, M., Hoekstra, M., Hildebrand, R.B., *et al.* *Arterioscler. Thromb. Vasc. Biol.* **27**, 2413-2419 (2007).
7. Doe, C., Bentley, R., Behm, D.J., *et al.* *J. Pharmacol. Exp. Ther.* **320**(1), 89-98 (2007).

Related Products

iPF_{2α}-IV - Item No. 16230 • 5-iPF_{2α}-VI - Item No. 16300 • iPF_{2α}-IV-d₄ - Item No. 316230 • iPF_{2α}-VI-d₄ - Item No. 316300 • 5-iPF_{2α}-VI-d₁₁ - Item No. 10006654 • 8,12-*iso*-iPF_{2α}-VI-d₁₁ - Item No. 10006878

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, 10/03/2011

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