

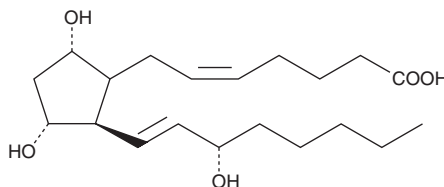
Product Information



Prostaglandin F_{2α}

Item No. 16010

CAS Registry No.: 551-11-1
Formal Name: 9α,11α,15S-trihydroxy-prosta-5Z,13E-dien-1-oic acid
Synonym: Dinoprost
MF: C₂₀H₃₄O₅
FW: 354.5
Purity: ≥99%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that Prostaglandin F_{2α} (PGF_{2α}) be stored as supplied at -20°C. It should be stable for at least two years.

PGF_{2α} is supplied as a crystalline solid. A stock solution may be made by dissolving the PGF_{2α} in an organic solvent purged with an inert gas. PGF_{2α} is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 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. Organic solvent-free aqueous solutions of PGF_{2α} can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of PGF_{2α} in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

PGF_{2α} is a widely distributed PG occurring in many species.¹⁻³ It causes contraction of vascular, bronchial, intestinal, and myometrial smooth muscle, and also exhibits potent luteolytic activity.¹ PGF_{2α} exhibits its receptor mediated physiological activity at 50-100 nM.¹ Maximal ovine myometrial contraction can be achieved at 125 nM PGF_{2α} *in vitro*.⁴

References

1. Samuelsson, B., Goldyne, M., Granström, E., *et al.* Prostaglandins and Thromboxanes. *Ann. Rev. Biochem.* **47**, 997-1029 (1978).
2. Speroff, L. and Ramwell, P.W. Prostaglandins in reproductive physiology. *Am. J. Obstet. Gynecol.* **107**, 1111-1130 (1970).
3. Watanabe, K., Iguchi, Y., Iguchi, S., *et al.* Stereospecific conversion of prostaglandin D₂ to (5Z,13E)-(15S)-9α,11β,15-trihydroxyprosta-5,13-dien-1-oic acid (9α,11β-prostaglandin F₂) and of prostaglandin H₂ to prostaglandin F_{2α} by bovine lung prostaglandin F synthase. *Proc. Natl. Acad. Sci. USA* **83**, 1583-1587 (1986).
4. Crankshaw, D.J. and Gaspar, V. Pharmacological characterization *in vitro* of prostanoid receptors in the myometrium of nonpregnant ewes. *J. Reprod. Fertil.* **103**, 55-61 (1995).

Related Products

Prostaglandin F_{2α} Alcohol - Item No. 16012 • Prostaglandin F_{2α} (tromethamine salt) - Item No. 16020 • Prostaglandin F_{2α} isopropyl ester - Item No. 16030 • Prostaglandin F_{2α} dimethyl amide - Item No. 16032 • Prostaglandin F_{2α} dimethyl amine - Item No. 16033 • Prostaglandin F_{2α}-d₄ - Item No. 316010 • Prostaglandin F_{2α} EIA Kit - Item No. 516011 • Prostaglandin F_{2α} EIA Kit (solid plate) - Item No. 516011.1 • COX (ovine) Inhibitor Screening Assay - Item No. 560101 • Chemiluminescent COX (ovine) Inhibitor Screening Assay - Item No. 760101 • Prostaglandin F_{2α} Lipid Maps MS Standard - Item No. 10007221 • (+)-Cloprostenol isopropyl ester - Item No. 10010016 • (+)-Cloprostenol methyl ester - Item No. 10010115 • (+)-Cloprostenol methyl amide - Item No. 10010495 • 17-phenoxy trinor Prostaglandin F_{2α} ethyl amide - Item No. 10010742

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, 04/08/2011

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