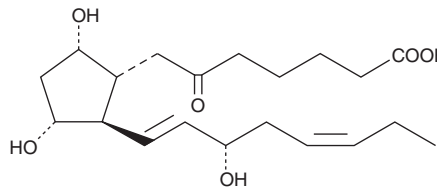


PRODUCT INFORMATION



Δ^{17} -6-keto Prostaglandin F_{1 α} Item No. 15230

CAS Registry No: 68324-95-8
Formal Name: 6-oxo-9 α ,11 α ,15S-trihydroxy-prosta-13E,17Z-dien-1-oic acid
Synonym: ω -3 6-keto PGF_{2 α}
MF: C₂₀H₃₂O₆
FW: 368.5
Purity: \geq 98%
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Δ^{17} -6-keto Prostaglandin F_{1 α} (Δ^{17} -6-keto PGF_{1 α}) 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 (DMF) purged with an inert gas can be used. The solubility of Δ^{17} -6-keto PGF_{1 α} in these is approximately >75 mg/ml and approximately >50 mg/ml in DMSO and DMF.

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 Δ^{17} -6-keto PGF_{1 α} is needed, the methyl acetate can be evaporated under a stream of nitrogen and the neat oil dissolved in the buffer of choice. Δ^{17} -6-keto PGF_{1 α} is soluble in PBS (pH 7.2) at a concentration of approximately 2.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Δ^{17} -6-keto PGF_{1 α} is a cyclooxygenase (COX) product of eicosapentaenoic acid (EPA) in various tissues such as seminal vesicles, lung, Polymorphonuclear leukocytes, and ocular tissues.^{1,2} Δ^{17} -6-keto PGF_{1 α} and other 3-series COX products from EPA, such as PGF_{3 α} , PGE₃, and thromboxane B₃, may be involved in the reduced incidence of glaucoma in patients on a marine-rich (EPA-rich) diet.²

References

1. Kulkarni, P.S., Kaufman, P.L., Srinivasan, B.D. Eicosapentaenoic acid metabolism in cynomolgus and rhesus conjunctiva and eyelid. *J. Ocul. Pharmacol.* **3**, 349-356 (1987).
2. Kulkarni, P.S. and Srinivasan, B.D. Eicosapentaenoic acid metabolism in human and rabbit anterior uvea. *Prostaglandins* **31**, 1159-1164 (1986).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/31/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM