

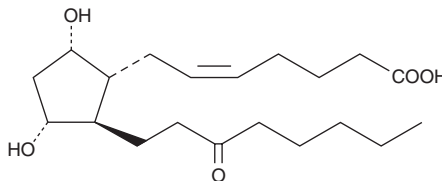
PRODUCT INFORMATION



13,14-dihydro-15-keto Prostaglandin F_{2α}

Item No. 16670

CAS Registry No.: 27376-76-7
Formal Name: 9α,11α-dihydroxy-15-oxo-prost-5-en-1-oic acid
Synonyms: 13,14-dihydro-15-keto PGF_{2α}, PGFM
MF: C₂₀H₃₄O₅
FW: 354.5
Purity: ≥95%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

13,14-dihydro-15-keto Prostaglandin F_{2α} (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 DMSO, dimethyl formamide, or ethanol purged with an inert gas can be used. The solubility of 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. Organic solvent-free solutions of 13,14-dihydro-15-keto PGF_{2α} can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 13,14-dihydro-15-keto 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.

Description

13,14-dihydro-15-keto PGF_{2α} is the first prominent plasma metabolite of PGF_{2α} in the 15-hydroxy PGDH pathway.¹ Measurement of 13,14-dihydro-15-keto PGF_{2α} in plasma can be used as a marker of the *in vivo* production of PGF_{2α}.¹⁻³

References

1. Samuelsson, B., Goldyne, M., Granström, E., *et al.* Prostaglandins and thromboxanes. *Annu. Rev. Biochem.* **47**, 997-1029 (1978).
2. Meyer, H.H.D., Eisele, K., and Osaso, J. A biotin-streptavidin amplified enzyme immunoassay for 13,14-dihydro-15-keto-PGF_{2α} Prostaglandins **38**, 375-383 (1989).
3. Del Vecchio, R.P., Maxey, K.M., and Lewis, G.S. A quantitative solid-phase enzyme immunoassay for 13,14-dihydro-15-keto-prostaglandin F_{2α} in plasma. *Prostaglandins* **43**, 321-330 (1992).

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, 01/19/2018

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