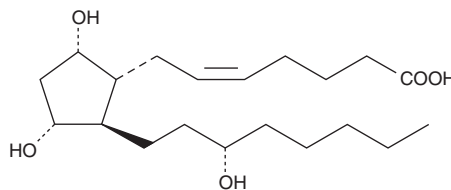


PRODUCT INFORMATION



13,14-dihydro Prostaglandin F_{2α} Item No. 16660

CAS Registry No.: 27376-74-5
Formal Name: 9α,11α,15S-trihydroxy-prost-5Z-en-1-oic acid
Synonym: 13,14-dihydro PGF_{2α}
MF: C₂₀H₃₆O₅
FW: 356.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that 13,14-dihydro prostaglandin F_{2α} (13,14-dihydro PGF_{2α}) be stored as supplied at -20°C. It should be stable for at least two years.

13,14-dihydro 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 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 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 PGF_{2α} in PBS (pH 7.2) is approximately 10 mg/ml. Store aqueous solutions of 13,14-dihydro PGF_{2α} on ice and use within 12 hours of preparation. Although the aqueous solutions of 13,14-dihydro PGF_{2α} may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Description

13,14-dihydro PGF_{2α} can cause luteolysis in hamsters with a potency five times that of PGF_{2α}. The ED₅₀ value for 13,14-dihydro PGF_{2α} as a luteolytic in hamsters is 1.5 μg/100 g¹

Reference

1. Andersen, N.H., Imamoto, S., Subramanian, N., *et al.* Molecular basis for prostaglandin potency. III. Tests of the significance of the "hairpin conformation" in biorecognition phenomena. *Prostaglandins* **22**, 841-856 (1981).

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, 08/13/2015

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