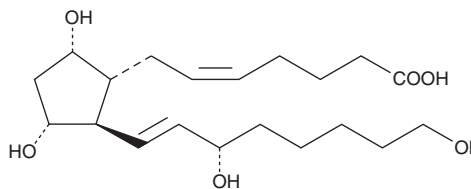


PRODUCT INFORMATION



20-hydroxy Prostaglandin F_{2α} Item No. 16950

CAS Registry No.: 57930-92-4
Formal Name: 9α,11α,15S,20-tetrahydroxy-prosta-5Z,13E-dien-1-oic acid
Synonym: 20-hydroxy PGF_{2α}
MF: C₂₀H₃₄O₆
FW: 370.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

20-hydroxy Prostaglandin F_{2α} (20-hydroxy 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, ethanol, and dimethyl formamide purged with an inert gas can be used. The solubility of 20-hydroxy 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. If an organic solvent-free solution of 20-hydroxy PGF_{2α} is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 20-hydroxy PGF_{2α} in PBS (pH 7.2) is approximately 10 mg/ml. Store aqueous solutions of 20-hydroxy PGF_{2α} on ice and use within 12 hours of preparation.

Description

20-hydroxy PGF_{2α} is the ω-oxidation product of PGF_{2α}. Cultured type II alveolar cells from pregnant rabbits metabolize exogenous PGF_{2α} via microsomal cytochrome P450 ω-oxidation, producing 20-hydroxy PGF_{2α} and its 15-hydroxy PGDH metabolites. Cells from male rabbits exhibit only the 15-hydroxy PGDH pathway.¹

Reference

1. Devereux, T.R., Fouts, J.R., and Eling, T.E. Metabolism of prostaglandin PG-F_{2α} by freshly isolated alveolar type II cells from lungs of adult male or pregnant rabbits. *Prostaglandins Leukot. Med.* **27**, 43-52 (1987).

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, 06/05/2019

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