

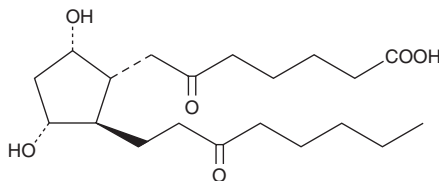
PRODUCT INFORMATION



6,15-diketo-13,14-dihydro Prostaglandin F_{1α}

Item No. 15270

CAS Registry No.: 63983-53-9
Formal Name: 6,15-dioxo-9α,11α-dihydroxy-prostan-1-oic acid
Synonym: 6,15-diketo-13,14-dihydro PGF_{1α}
MF: C₂₀H₃₄O₆
FW: 370.5
Purity: ≥98% (mixture of tautomers)
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

6,15-diketo-13,14-dihydro PGF_{1α} 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, or dimethyl formamide purged with an inert gas can be used. The solubility of 6,15-diketo-13,14-dihydro PGF_{1α} in these solvents is approximately 16, 50, and 100 mg/ml, respectively.

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 aqueous solution of 6,15-diketo-13,14-dihydro PGF_{1α} is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 6,15-diketo-13,14-dihydro PGF_{1α} in PBS (pH 7.2) is approximately 3.3 mg/ml. Store aqueous solutions of 6,15-diketo-13,14-dihydro PGF_{1α} on ice and use within 12 hours of preparation.

Description

6,15-diketo-13,14-dihydro PGF_{1α} is a metabolite of PGI₂. It was shown to enhance intracellular cAMP and cholesterol catabolism in bovine arterial smooth muscle cells.¹

Reference

1. Etingin, O.R., Weksler, B.B., and Hajjar, D.P. Cholesterol metabolism is altered by hydrolytic metabolites of prostacyclin in arterial smooth muscle cells. *J. Lipid Res.* **27**, 530-536 (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, 03/14/2024

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