

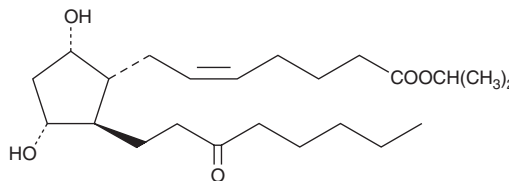
PRODUCT INFORMATION



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

Item No. 16671

CAS Registry No.: 199920-18-8
Formal Name: 9α,11α-dihydroxy-15-oxo-Prost-5Z-en-1-oic acid, 1-methylethyl ester
Synonym: 13,14-dihydro-15-keto PGF_{2α} isopropyl ester
MF: C₂₃H₄₀O₅
FW: 396.6
Purity: ≥98%
Supplied as: A 1 mg/ml 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

13,14-dihydro-15-keto Prostaglandin F_{2α} isopropyl ester (13,14-dihydro-15-keto PGF_{2α} isopropyl ester) 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 13,14-dihydro-15-keto PGF_{2α} isopropyl ester in ethanol and DMSO is approximately 50 mg/ml and approximately 30 mg/ml in DMF.

13,14-dihydro-15-keto PGF_{2α} isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 13,14-dihydro-15-keto PGF_{2α} isopropyl ester should be diluted with the aqueous buffer of choice. The solubility of 13,14-dihydro-15-keto PGF_{2α} isopropyl ester in PBS (pH 7.2) is approximately 0.05 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

13,14-dihydro-15-keto Prostaglandin F_{2α} (PGF_{2α}) isopropyl ester is an esterified form of 13,14-dihydro-15-keto PGF_{2α} (Item No. 16670).

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, 12/18/2025

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