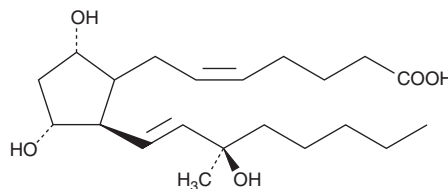


PRODUCT INFORMATION



15(R)-15-methyl Prostaglandin F_{2α} Item No. 16730

CAS Registry No.: 35864-81-4
Formal Name: 9α,11α,15R-trihydroxy-15-methyl-prosta-5Z,13E-dien-1-oic acid
Synonym: 15(R)-15-methyl PGF_{2α}
MF: C₂₁H₃₆O₅
FW: 368.5
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

15(R)-15-methyl Prostaglandin F_{2α} (15(R)-15-methyl 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 15(R)-15-methyl PGF_{2α} in these solvents is approximately 100 mg/ml. The solubility of 15(R)-15-methyl PGF_{2α} in 10 mM Na₂CO₃ is approximately 6.5 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 15(R)-15-methyl 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 15(R)-15-methyl 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

15(R)-15-methyl PGF_{2α} is a metabolically stable analog of PGF_{2α}. 15(R)-15-methyl PGF_{2α} is an inactive, prodrug PGF agonist designed for activation by gastric acid after oral administration. Acid-catalyzed epimerization of 15(R)-15-methyl PGF_{2α} converts it into the active 15(S)-isomer.^{1,2} The 15(S)-isomer induces luteolysis when injected in rhesus monkeys at a dose of about 12 mg/animal, while the 15(R)-isomer does not.³

References

1. Plaisted, S.M., DeZwaan, J., and Snider, B.G. High-performance liquid chromatographic determination of acid-catalyzed degradation products of methyl carboprost in a polymeric controlled-release device. *J. Chromatogr.* **314**, 369-377 (1984).
2. Hamberg, M., Zhang, L.Y., and Bergström, S. On the pH-dependent degradation of 15(S)-15 methyl-prostaglandin F_{2α} (Carboprost). *Eur. J. Pharm. Sci.* **3(1)**, 27-38 (1995).
3. Wilks, J.W. Inhibition of the monkey corpus luteum with 15-methyl prostaglandins. *Prostaglandins* **20(5)**, 793-805 (1980).

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, 03/25/2024

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