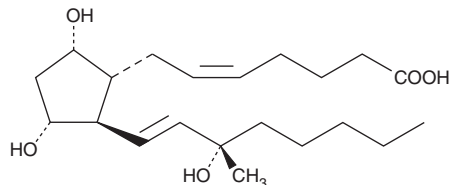


PRODUCT INFORMATION



15(S)-15-methyl Prostaglandin F_{2α} Item No. 16743

CAS Registry No.: 35700-23-3
Formal Name: 9α,11α,15S-trihydroxy-15-methyl-prosta-5Z,13E-dien-1-oic acid
Synonyms: Carboprost, 15(S)-15-methyl PGF_{2α}, U-32921
MF: C₂₁H₃₆O₅
FW: 368.5
Purity: ≥98%
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(S)-15-methyl Prostaglandin F_{2α} (15(S)-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(S)-15-methyl PGF_{2α} in these solvents is approximately 100 mg/ml.

15(S)-15-methyl PGF_{2α} is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15(S)-15-methyl PGF_{2α} should be diluted with the aqueous buffer of choice. The solubility of 15(S)-15-methyl PGF_{2α} in PBS (pH 7.2) and 10 mM Na₂CO₃ is approximately 10 and 6.5 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

15(S)-15-methyl PGF_{2α} is a metabolically stable analog of PGF_{2α}. It is a potent uterine stimulant and abortifacient which can be administered intramuscularly to induce labor.¹⁻³ It induces luteolysis and reduces serum progesterone concentrations when given as an intramuscular injection in non-pregnant rhesus monkeys.⁴

References

1. Carlan, S.J., Gushwa, J.P., O'Brien, W.F., *et al.* Effect of intramuscular 15-methyl prostaglandin F_{2α} after second-trimester delivery. *Obstet. Gynecol.* **89(1)**, 5-9 (1997).
2. Lindblom, B., Källfelt, B., Hahlin, M., *et al.* Local prostaglandin F_{2α} injection for termination of ectopic pregnancy. *Lancet* **1(8536)**, 776-777 (1987).
3. Adaikan, P.G., Prasad, R.N.V., Kottegoda, S.R., *et al.* Effect of indomethacin on the injection-abortion interval of 15(S)15 methyl PGF_{2α}-induced mid-trimester abortions - a randomized study. *Prostaglandins Leukot. Med.* **27(2-3)**, 161-167 (1987).
4. 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, 03/15/2024

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