

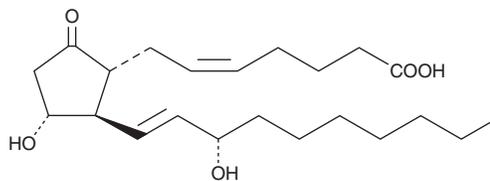
PRODUCT INFORMATION



20-ethyl Prostaglandin E₂

Item No. 14940

CAS Registry No.: 37492-24-3
Formal Name: (5Z)-7-[(1R,2R,3R)-3-hydroxy-2-[(1E,3S)-3-hydroxy-1-decen-1-yl]-5-oxocyclopentyl]-5-heptenoic acid
Synonym: 20-ethyl PGE₂
MF: C₂₂H₃₆O₅
FW: 380.5
Purity: ≥97%
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

20-ethyl Prostaglandin E₂ (20-ethyl PGE₂) 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 20-ethyl PGE₂ in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

20-ethyl PGE₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 20-ethyl PGE₂ should be diluted with the aqueous buffer of choice. The solubility of 20-ethyl PGE₂ in PBS (pH 7.2) is approximately 0.125 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

20-ethyl PGE₂ is an analog of PGE₂ in which the ω-chain has been extended by the addition of two methylene carbon atoms. The only well studied prostaglandin analog with this structural feature is unoprostone, an F-series prostaglandin that is clinically approved as a glaucoma medication.¹ Unoprostone also contains lower side chain modifications (13,14-dihydro-15-keto) which severely limit its affinity for FP receptors, contributing to its lack of potency as a medication. 20-ethyl PGE₂ retains the natural 15(S) allylic hydroxyl in the lower side chain, which may improve its potency relative to unoprostone. However, ligand binding assays of this analog with respect to EP or other prostanoid receptors have not been published. E-type prostaglandins have been widely reported to have inflammatory, cytoprotective, and a variety of other effects.²⁻⁴

References

1. Hara, M. and Spencer, C.M. Unoprostone (isopropyl unoprostone). *Drugs Aging* **9**(3), 213-218 (1996).
2. Matsumoto, H., Naraba, H., Murakami, M., *et al.* Concordant induction of prostaglandin E₂ synthase with cyclooxygenase-2 leads to preferred production of prostaglandin E₂ over thromboxane and prostaglandin D₂ in lipopolysaccharide-stimulated rat peritoneal macrophages. *Biochem. Biophys. Res. Commun.* **230**(1), 110-114 (1997).
3. Karim, S.M.M., Carter, D.C., Bhana, D., *et al.* Effect of orally administered prostaglandin E₂ and its 15-methyl analogues on gastric secretion. *Br. Med. J.* **1**(5846), 143-146 (1973).
4. Jackson, G.M., Sharp, H.T., and Varner, M.W. Cervical ripening before induction of labor: A randomized trial of prostaglandin E₂ gel versus low-dose oxytocin. *Am. J. Obstet. Gynecol.* **171**(4), 1092-1096 (1994).

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, 03/13/2024

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