

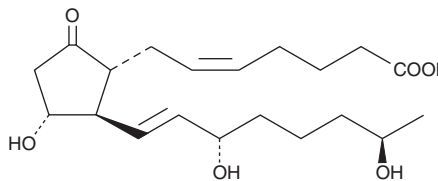
PRODUCT INFORMATION



19(R)-hydroxy Prostaglandin E₂

Item No. 14910

CAS Registry No.: 64625-54-3
Formal Name: 11 α ,15S,19R-trihydroxy-9-oxo-prosta-5Z,13E-dien-1-oic acid
Synonyms: 19(R)-hydroxy PGE₂, 19(R)-OH PGE₂
MF: C₂₀H₃₂O₆
FW: 368.5
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

19(R)-hydroxy Prostaglandin E₂ (19(R)-hydroxy PGE₂) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol 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 19(R)-hydroxy PGE₂ in these solvents is approximately 100 mg/ml.

19(R)-hydroxy PGE₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 19(R)-hydroxy PGE₂ should be diluted with the aqueous buffer of choice. The solubility of 19(R)-hydroxy PGE₂ in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

19(R)-hydroxy PGE₂ is found in the semen of primates, including man.¹ It is a potent smooth muscle relaxant and a selective agonist for the EP₂ receptor.^{1,2} It has an EC₅₀ value of 200 nM for relaxing cat tracheal rings which express EP₂ receptors.²

References

1. Kelly, R.W., Taylor, P.L., Hearn, J.P., *et al.* 19-Hydroxyprostaglandin E₁ as a major component of the semen of primates. *Nature* **260(5551)**, 544-545 (1976).
2. Woodward, D.F., Protzman, C.E., Krauss, A.H.P., *et al.* Identification of 19(R)-OH prostaglandin E₂ as a selective prostanoid EP₂-receptor agonist. *Prostaglandins* **46(4)**, 371-383 (1993).

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, 02/06/2024

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