

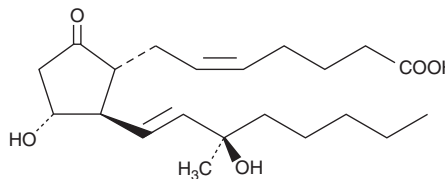
PRODUCT INFORMATION



15(R)-15-methyl Prostaglandin E₂

Item No. 14725

CAS Registry No.: 55028-70-1
Formal Name: 9-oxo-11 α ,15R-dihydroxy-15-methyl-prosta-5Z,13E-dien-1-oic acid
Synonyms: Arbaprostil, 15(R)-15-methyl PGE₂
MF: C₂₁H₃₄O₅
FW: 366.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

15(R)-15-methyl Prostaglandin E₂ (15(R)-15-methyl PGE₂) is supplied as a crystalline solid. A stock solution may be made by dissolving the 15(R)-15-methyl PGE₂ in the solvent of choice, which should be purged with an inert gas. 15(R)-15-methyl PGE₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 15(R)-15-methyl PGE₂ in these solvents is approximately 100 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. Organic solvent-free aqueous solutions of 15(R)-15-methyl PGE₂ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 15(R)-15-methyl 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

15(R)-15-methyl PGE₂ is a prodrug for the potent PGE₂ (Item No. 14010) analog 15(S)-15-methyl PGE₂ (Item No. 14730).¹ Acid-catalyzed epimerization in the stomach produces the 15(S)-hydroxy compound which is biologically active.² Oral administration of 15(R)-15-methyl PGE₂ to dogs or rats at 10-300 μ g/kg results in a dose-dependent inhibition of gastric acid secretion and an increase in the rate of duodenal bicarbonate secretion.^{3,4}

References

1. Yankee, E.W., Axen, U., and Bundy, G.L. Total synthesis of 15-methylprostaglandins. *J. Am. Chem. Soc.* **96**, 5865-5876 (1974).
2. Takanashi, H., Kawabe, Y., and Akima, M. Acid-promoted epimerization of arbaprostil, 15(R)-15-methylprostaglandin E₂, elicits gastric antisecretory activities in rats. *Jpn. J. Pharmacol.* **57**, 559-564 (1991).
3. Takanashi, H. and Itoh, Z. Gastric antisecretory activity of 15(R)-15-methylprostaglandin E₂, arbaprostil, in dogs. *Jpn. J. Pharmacol.* **57**, 447-451 (1991).
4. Li, J., Nagata, T., Yoshida, M., et al. Effect of 15(R)-15-methyl PGE₂ (arbaprostil) on duodenal bicarbonate secretion in rat. *Gastroenterol. Jpn.* **24**, 8-11 (1989).

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, 01/31/2024

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