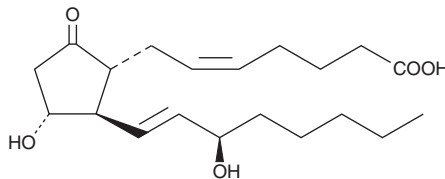


PRODUCT INFORMATION

15(R)-Prostaglandin E₂

Item No. 14710

CAS Registry No.: 38873-82-4
Formal Name: 11 α ,15R-dihydroxy-9-oxo-prosta-5Z,13E-dien-1-oic acid
Synonyms: 15-*epi* PGE₂, 15(R)-PGE₂
MF: C₂₀H₃₂O₅
FW: 352.5
Purity: $\geq 98\%$
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(R)-Prostaglandin E₂ (15(R)-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 purged with an inert gas can be used. The solubility of 15(R)-PGE₂ in these solvents is approximately 100 mg/ml.

15(R)-PGE₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15(R)-PGE₂ should be diluted with the aqueous buffer of choice. The solubility of 15(R)-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)-PGE₂ is the opposite C-15 stereoisomer of the 15(S)-PGE₂. 15(R)-PGE₂ is much less potent than the natural stereoisomer in most biological assays. In hamsters, a dose of 0.20 mg of PGE₂ (Item No. 14010) inhibits fertility by 100%, whereas a dose of 1.0 mg of 15(R)-PGE₂ inhibits fertility by only 17%.¹ 15(R)-PGE₂ occurs as a minor impurity in some pharmaceutical formulations of PGE₂ used in obstetrics.

Reference

1. Miller, W.L. and Sutton, M.J. Relative biological activity of certain prostaglandins and their enantiomers. *Prostaglandins* **11**(1), 77-84 (1976).

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