

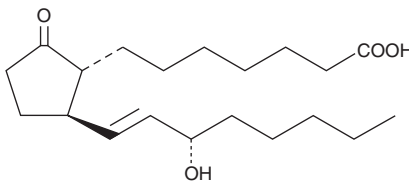
PRODUCT INFORMATION



11-deoxy Prostaglandin E₁

Item No. 13510

CAS Registry No.: 37786-00-8
Formal Name: 9-oxo-15S-hydroxy-prost-13E-en-1-oic acid
Synonym: 11-deoxy PGE₁
MF: C₂₀H₃₄O₄
FW: 338.5
Purity: ≥96%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

11-deoxy Prostaglandin E₁ (11-deoxy PGE₁) is supplied as a crystalline solid. A stock solution may be made by dissolving the 11-deoxy PGE₁ in the solvent of choice, which should be purged with an inert gas. 11-deoxy PGE₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 11-deoxy PGE₁ in ethanol and DMSO is approximately 50 mg/ml and approximately 100 mg/ml in DMF.

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 11-deoxy PGE₁ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 11-deoxy PGE₁ in PBS (pH 7.2) is approximately 1.6 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

11-deoxy PGE₁ is a synthetic analog of PGE₁. Early reports show that it is a selective agonist for the EP₂ receptor but effective at much higher concentrations than PGE₂.^{1,2} However, later studies show that it is a non-selective agonist of EP receptors and stimulates cAMP release in Jurkat cells with an EC₅₀ of 0.25 μM.³ 11-deoxy PGE₁ also exhibits vasodepressor and bronchodilator responses in guinea pigs.⁴ 11-deoxy PGE₁ exhibits selectivity for the mouse EP₃ receptor and is essentially equipotent to PGE₁ at this receptor subtype. The K_i values for binding to the mouse EP₁, EP₂, EP₃, and EP₄ receptors are 600, 45, 1.1, and 23 nM, respectively.⁵

References

1. Dong, Y.J., Jones, R.L. and Wilson, N.H. Prostaglandin E receptor subtypes in smooth muscle: Agonist activities of stable prostacyclin analogues. *Br. J. Pharmacol.* **87(1)**, 97-107 (1986).
2. Chen, J. and Woodward, D.F. Prostanoid-induced relaxation of precontracted cat ciliary muscle is mediated by EP₂ and DP receptors. *Invest. Ophthalmol. Vis. Sci.* **33(11)**, 3195-3201 (1992).
3. De Vries, G.W., Guarino, P., McLaughlin, A., et al. An EP receptor with a novel pharmacological profile in the T-cell line Jurkat. *Br. J. Pharmacol.* **115(7)**, 1231-1234 (1995).
4. Hall, D.W.R. and Jaitly, K.D. Structure-activity relationships in a series of 11-deoxy prostaglandins. *Prostaglandins* **11(3)**, 573-587 (1976).
5. Kiriya, M., Ushikubi, F., Kobayashi, T., et al. Ligand binding specificities of the eight types and subtypes of the mouse prostanoid receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **122(2)**, 217-224 (1997).

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, 03/21/2024

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