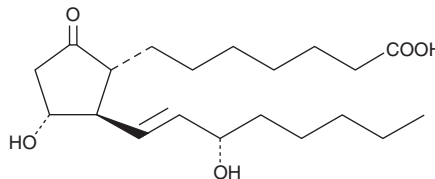


PRODUCT INFORMATION



Prostaglandin E₁ Item No. 13010

CAS Registry No.: 745-65-3
Formal Name: 11 α ,15S-dihydroxy-9-oxo-prost-13E-en-1-oic acid
Synonyms: Alprostadil, NSC 165559, PGE₁
MF: C₂₀H₃₄O₅
FW: 354.5
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

Prostaglandin E₁ (PGE₁) is supplied as a crystalline solid. A stock solution may be made by dissolving the PGE₁ in the solvent of choice, which should be purged with an inert gas. PGE₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 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 PGE₁ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of PGE₁ in PBS (pH 7.2) is approximately 0.5 mg/ml. Avoid adding PGE₁ to basic solutions (pH >7.4), since base treatment will degrade PGE₁ to PGA and PGB compounds. We do not recommend storing the aqueous solution for more than one day.

Description

PGE₁ is a vasoactive prostaglandin and an active metabolite of dihomo- γ -linolenic acid (DGLA; Item No. 90230).^{1,2} It is formed from DGLA by COX-1 and COX-2. PGE₁ is an agonist of the PGE₂ (Item No. 14010) receptor subtypes EP₁, EP₂, EP₃, and EP₄, and the IP receptor (K_s = 36, 10, 1.1, 2.1, and 33 nM, respectively, for the mouse receptors).³ It inhibits ADP-induced platelet aggregation of isolated human platelet-rich plasma (IC₅₀ = 40 nM) and isoproterenol-induced increases in L-type calcium current (I_{Ca}) in isolated rabbit atrial cells (EC₅₀ = 27 nM).^{4,5} PGE₁ (100 nM) induces vasodilation in isolated rat aortic rings and activates ATP-sensitive potassium channels (K_{ATP}) in a cell-attached patch clamp assay using isolated rat vascular smooth muscle cells (VSMCs).¹ It decreases femoral arterial perfusion pressure in dogs.⁶ Formulations containing PGE₁ have been used in the treatment of erectile dysfunction and to maintain patency of the ductus arteriosus in neonates with congenital heart defects who depend on a patent ductus arteriosus for survival.

References

1. Eguchi, S., Kawano, T., Yinhu, *et al.* *J. Cardiovasc. Pharmacol.* **50(6)**, 686-691 (2007).
2. Levin, G., Duffin, K.L., Obukowicz, M.G., *et al.* *Biochem. J.* **365(Pt 2)**, 489-496 (2002).
3. Kiriya, M., Ushikubi, F., Kobayashi, T., *et al.* *Br. J. Pharmacol.* **122(2)**, 217-224 (1997).
4. Kobzar, G., Mardla, V., Järving, I., *et al.* *Proc. Estonian Acad. Sci. Chem.* **40(N3)**, 179-180 (1991).
5. Yamamoto, T., Habuchi, Y., Tanaka, H., *et al.* *Am. J. Physiol.* **277(4)**, H1369-H1374 (1999).
6. Nakano, J. *Br. J. Pharmacol.* **44(1)**, 63-70 (1972).

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, 04/12/2023

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