

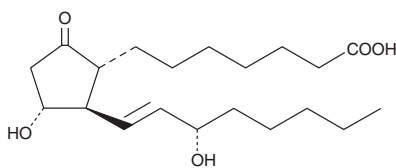
PRODUCT INFORMATION



Prostaglandin E₁ Quant-PAK Item No. 10006844

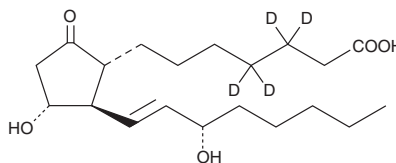
Prostaglandin E₁

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%
Stability: \geq 4 years at -20°C
Supplied as: A crystalline solid



Prostaglandin E₁-d₄

CAS Registry No.: 211105-33-8
Formal Name: 9-oxo-11 α ,15S-dihydroxy-prost-13E-en-1-oic-3,3,4,4-d₄ acid
Synonyms: Alprostadil-d₄, PGE₁-d₄
MF: C₂₀H₃₀D₄O₅
FW: 358.5
Chemical Purity: \geq 98% (Prostaglandin E₁)
Deuterium Incorporation: \geq 99% deuterated forms (d₁-d₄); \leq 1% d₀
Stability: \geq 4 years at -20°C
Supplied as: A solution in methyl acetate



This prostaglandin E₁ (PGE₁) Quant-PAK contains 50 μ g of PGE₁-d₄ and 2-4 mg of PGE₁ (please see the vial for exact amount and concentration).

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.

PGE₁-d₄ 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 (DMF) purged with an inert gas can be used. The solubility of PGE₁-d₄ in ethanol and DMSO is approximately 50 mg/ml and approximately 100 mg/ml in DMF.

PGE₁-d₄ contains four deuterium atoms at the 3, 3', 4, and 4' positions. It is intended for use as an internal standard for the quantification of PGE₁ by GC- or LC-mass spectrometry. The accuracy of the sample weight in the PGE₁-d₄ vial is between 5% over and 2% under the weight indicated on the vial. For better precision we have provided a precisely weighed unlabeled PGE₁, with the precise weight (2-4 mg) indicated on the vial. Using this vial the deuterated standard can be quantified by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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_is = 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, 11/10/2023

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