

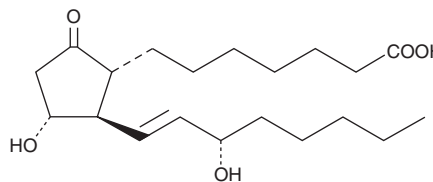
# Product Information



## Prostaglandin E<sub>1</sub>

Item No. 13010

**CAS Registry No.:** 745-65-3  
**Formal Name:** 9-oxo-11 $\alpha$ ,15S-dihydroxy-prost-13E-en-1-oic acid  
**Synonym:** Alprostadil  
**MF:** C<sub>20</sub>H<sub>34</sub>O<sub>5</sub>  
**FW:** 354.5  
**Purity:**  $\geq$ 98%  
**Stability:**  $\geq$ 2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

Prostaglandin E<sub>1</sub> (PGE<sub>1</sub>) is a clinically useful vasodilator and an inhibitor of platelet aggregation. For long term storage, we suggest that PGE<sub>1</sub> be stored as supplied at -20°C. It should be stable for at least two years.

PGE<sub>1</sub> is supplied as a crystalline solid. A stock solution may be made by dissolving the PGE<sub>1</sub> in an organic solvent. PGE<sub>1</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PGE<sub>1</sub> in these solvents is approximately 50 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 PGE<sub>1</sub> can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of PGE<sub>1</sub> in PBS (pH 7.2) is approximately 1.67 mg/ml. Avoid adding PGE<sub>1</sub> to basic solutions (pH > 7.4), since base treatment will degrade PGE<sub>1</sub> to PGA and PGB compounds. We do not recommend storing the aqueous solution for more than one day.

PGE<sub>1</sub> is the theoretical cyclooxygenase metabolite of dihomo- $\gamma$ -linolenic acid (DGLA), but it is virtually undetectable in the plasma of normal humans or other animals.<sup>1</sup> Its pharmacology includes vasodilation, hypotension, and anti-platelet activities. The IC<sub>50</sub> of PGE<sub>1</sub> for the inhibition of ADP-induced human platelet aggregation is 40 nM.<sup>2,3</sup> The vasorelaxant and anti-hypertensive effects of PGE<sub>1</sub> are used to treat male erectile dysfunction and to provide emergency vasodilation of the patent ductus arteriosus in infants whose cardiac anomalies require pulmonary shunting for survival.<sup>4,5</sup> In human males, the intracavernosal effective dose range for PGE<sub>1</sub> is 2 to 80  $\mu$ g, and the transurethral range is 125 to 1,000  $\mu$ g.<sup>4</sup>

### References

1. Cawello, W., Schweer, H., Dietrich, B., *et al.* Pharmacokinetics of prostaglandin E<sub>1</sub> and its main metabolites after intracavernous injection and short-term infusion of prostaglandin E<sub>1</sub> in patients with erectile dysfunction. *J. Urol.* **158**, 1403-1407 (1997).
2. Kobzar, G., Mardla, V., Järving, I., *et al.* Antiaggregating potency of E-type prostaglandins in human and rabbit platelets. *Proc. Estonian Acad. Sci. Chem.* **40**, 179-180 (1991).
3. Okada, F., Nukada, T., Yamauchi, Y., *et al.* The hypotensive effect of prostaglandin E<sub>1</sub> on hypertensive cases of various types. *Prostaglandins* **7**, 99-106 (1974).
4. Padma-Nathan, H., Hellstrom, W.J.G., Kaiser, F.E., *et al.* Treatment of men with erectile dysfunction with transurethral alprostadil. *N. Engl. J. Med.* **336**, 1-7 (1997).
5. Olley, P.M. and Cocceani, F. Prostaglandins and the ductus arteriosus. *Annu. Rev. Med.* **32**, 375-3785 (1981).

### Related Products

Prostaglandin E<sub>1</sub> (analytical grade)- Item No. 13010.1 • Prostaglandin E<sub>1</sub> (GMP,USP)- Item No. 13010.5

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

**For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 06/20/2011

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com