

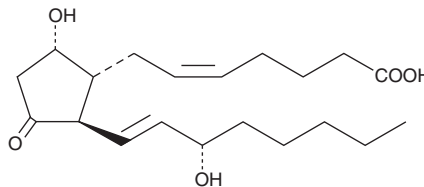
# PRODUCT INFORMATION



## Prostaglandin D<sub>2</sub>

Item No. 12010

CAS Registry No.: 41598-07-6  
Formal Name: 9 $\alpha$ ,15S-dihydroxy-11-oxo-prosta-5Z,13E-dien-1-oic acid  
Synonym: PGD<sub>2</sub>  
MF: C<sub>20</sub>H<sub>32</sub>O<sub>5</sub>  
FW: 352.5  
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 D<sub>2</sub> (PGD<sub>2</sub>) is supplied as a crystalline solid. A stock solution may be made by dissolving the PGD<sub>2</sub> in the solvent of choice, which should be purged with an inert gas. PGD<sub>2</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PGD<sub>2</sub> in these solvents is approximately 75, 50, and 100 mg/ml, respectively.

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 PGD<sub>2</sub> can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of PGD<sub>2</sub> in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PGD<sub>2</sub> is the major eicosanoid product of mast cells and is released in large quantities during allergic and asthmatic anaphylaxis.<sup>1</sup> Mastocytosis patients produce excessive amounts of PGD<sub>2</sub>, which causes vasodilation, flushing, hypotension, and syncopal episodes.<sup>1</sup> PGD<sub>2</sub> is also produced in the brain via an alternative pathway involving a soluble, secreted PGD-synthase also known as  $\beta$ -trace.<sup>2,3</sup> In the brain, PGD<sub>2</sub> produces normal physiological sleep and lowering of body temperature.<sup>2,3</sup> Further pharmacological actions include inhibition of platelet aggregation and relaxation of vascular smooth muscle.<sup>4</sup> PGD<sub>2</sub> inhibits human ovarian tumor cell proliferation with an IC<sub>50</sub> of 6.8  $\mu$ M.<sup>5</sup>

### References

1. Roberts, L.J., II, and Sweetman, B.J. Metabolic fate of endogenously synthesized prostaglandin D<sub>2</sub> in a human female with mastocytosis. *Prostaglandins* **30(3)**, 383-400 (1985).
2. Hayaishi, O. Sleep-wake regulation by prostaglandins D<sub>2</sub> and E<sub>2</sub>. *J. Biol. Chem.* **263(29)**, 14593-14596 (1988).
3. Onoe, H., Ueno, R., Fujita, I., et al. Prostaglandin D<sub>2</sub>, a cerebral sleep-inducing substance in monkeys. *Proc. Natl. Acad. Sci. USA* **85(11)**, 4082-4086 (1988).
4. Giles, H. and Leff, P. The biology and pharmacology of PGD<sub>2</sub>. *Prostaglandins* **35(2)**, 277-300 (1988).
5. Kikuchi, Y., Kita, T., Hirata, J., et al. Preclinical studies of antitumor prostaglandins by using human ovarian cancer cells. *Cancer Metastasis Rev.* **13(3-4)**, 309-315 (1994).

#### 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, 02/26/2024

#### 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