

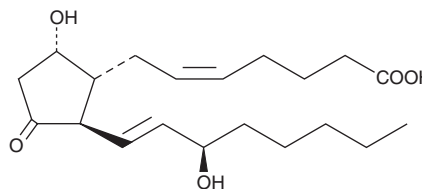
PRODUCT INFORMATION



15(R)-Prostaglandin D₂

Item No. 10118

CAS Registry No.: 59894-05-2
Formal Name: 9 α ,15R-dihydroxy-11-oxo-prosta-5Z,13E-dien-1-oic acid
Synonym: 15(R)-PGD₂
MF: C₂₀H₃₂O₅
FW: 352.5
Purity: \geq 98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

15(R)-Prostaglandin D₂ (15(R)-PGD₂) 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 purged with an inert gas can be used. The solubility of 15(R)-PGD₂ in these solvents is approximately 50 mg/ml.

15(R)-PGD₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15(R)-PGD₂ should be diluted with the aqueous buffer of choice. The solubility of 15(R)-PGD₂ in PBS (pH 7.4) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Many of the effects of PGD₂ are transduced *via* a traditional 7-transmembrane GPCR, the DP₁ receptor.¹ However, in certain leukocytes and other immune cells, a second PGD₂ receptor referred to as the CRTH2 or DP₂ receptor has been cloned and characterized.^{2,3} 15(R)-15-methyl PGD₂ is a suprisingly potent agonist at the DP₂ receptor, being about five times more potent than PGD₂ itself.⁴ 15(R)-PGD₂ has the potential to be produced endogenously as a natural ligand for DP₂, unlike the synthetic 15-methyl analogs. In preliminary reports, 15(R)-PGD₂ has also been reported to be a potent and selective DP₂ receptor agonist.^{5,6}

References

1. Boie, Y., Sawyer, N., Slipetz, D.M., *et al.* *J. Biol. Chem.* **270**(32), 18910-18916 (1995).
2. Abe, H., Takeshita, T., Nagata, K., *et al.* *Gene* **227**(1), 71-77 (1999).
3. Hirai, H., Tanaka, K., Yoshie, O., *et al.* *J. Exp. Med.* **193**(2), 255-261 (2001).
4. Monneret, G., Cossette, C., Gravel, S., *et al.* *J. Pharmacol. Exp. Ther.* **304**(1), 349-355 (2003).
5. Kim, S., Bellone, S., Maxey, K.M., *et al.* *Bioorganic & Medicinal Chemistry Letters* **15**(7), 1873-1876 (2005).
6. Cossette, C., Walsh, S.E., Kim, S., *et al.* *J. Pharmacol. Exp. Ther.* **320**(1), 173-179 (2007).

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, 02/27/2024

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