

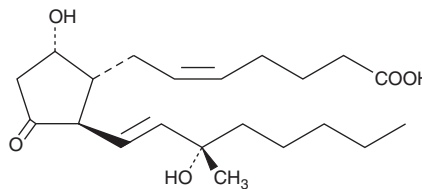
PRODUCT INFORMATION



15(S)-15-methyl Prostaglandin D₂

Item No. 12730

CAS Registry No.: 85280-90-6
Formal Name: 9 α ,15S-dihydroxy-11-oxo-15-methyl-prosta-5Z,13E-dien-1-oic acid
Synonym: 15(S)-15-methyl PGD₂
MF: C₂₁H₃₄O₅
FW: 366.5
Purity: ≥95%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

15(S)-15-methyl Prostaglandin D₂ (15(S)-15-methyl 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(S)-15-methyl PGD₂ in these solvents is approximately 75, 50, and 100 mg/ml, respectively.

15(S)-15-methyl PGD₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15(S)-15-methyl PGD₂ should be diluted with the aqueous buffer of choice. The solubility of 15(S)-15-methyl PGD₂ in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15(S)-15-methyl PGD₂ is a metabolically stable synthetic analog of PGD₂ (Item No. 12010). In contrast to PGD₂, 15(S)-15-methyl PGD₂ induces vasoconstriction and increases systemic blood pressure with much reduced inhibitory activity on ADP-induced platelet aggregation.¹ It also exhibits strong antifertility activity in hamsters (200-fold more potent than PGD₂).¹

Reference

1. Bundy, G.L., Morton, D.R., Peterson, D.C., *et al.* Synthesis and platelet aggregation inhibiting activity of prostaglandin D analogues. *J. Med. Chem.* **26**(6), 790-799 (1983).

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, 03/08/2024

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