

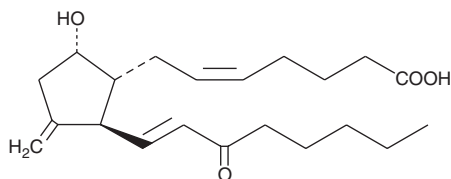
PRODUCT INFORMATION



11-deoxy-11-methylene-15-keto Prostaglandin D₂

Item No. 12415

CAS Registry No.: 958759-75-6
Formal Name: 9 α -hydroxy-11-methylene-15-oxo-prosta-5Z,13E-dien-1-oic acid
Synonym: 11-deoxy-11-methylene-15-keto PGD₂
MF: C₂₁H₃₂O₄
FW: 348.5
Purity: $\geq 98\%$
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

11-deoxy-11-methylene-15-keto Prostaglandin D₂ 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 11-deoxy-11-methylene-15-keto prostaglandin D₂ in these solvents is approximately 100 mg.

11-deoxy-11-methylene-15-keto Prostaglandin D₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the aqueous solution of 11-deoxy-11-methylene-15-keto prostaglandin D₂ should be diluted with the aqueous buffer of choice. The solubility of 11-deoxy-11-methylene-15-keto prostaglandin D₂ in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

11-deoxy-11-methylene-15-keto Prostaglandin D₂ is a stabilized form of 15-keto prostaglandin D₂ (15-keto PGD₂).

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, 07/21/2026

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