

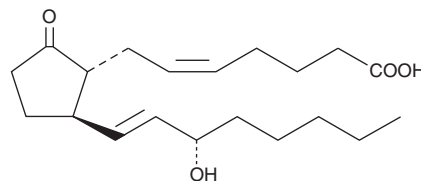
PRODUCT INFORMATION



11-deoxy Prostaglandin E₂

Item No. 14520

CAS Registry No.: 35536-53-9
Formal Name: 15S-hydroxy-9-oxo-prosta-5Z,13E-dien-1-oic acid
Synonym: 11-deoxy PGE₂
MF: C₂₀H₃₂O₄
FW: 336.5
Purity: ≥99%
Supplied as: A crystalline solid
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 Prostaglandin E₂ (11-deoxy PGE₂) is supplied as a crystalline solid. A stock solution may be made by dissolving the 11-deoxy PGE₂ in the solvent of choice, which should be purged with an inert gas. 11-deoxy PGE₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 11-deoxy PGE₂ in these solvents is approximately 100 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 11-deoxy PGE₂ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 11-deoxy PGE₂ in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

11-deoxy PGE₂ is a stable, synthetic analog of PGE₂ (Item No. 14010). In contrast to PGE₂ which has bronchodilation effects, 11-deoxy PGE₂ is a powerful bronchoconstrictor and contracts human respiratory tract smooth muscle with potencies ranging from 5 to 30 times higher than PGF_{2α} (Item No. 16010).¹

Reference

1. Karim, S.M.M., Adaikan, P.G., and Kottogoda, S.R. Prostaglandins and human respiratory tract smooth muscle: Structure activity relationship. *Adv. Prostaglandin Thromboxane Res.* **7**, 969-980 (1980).

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, 04/02/2024

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