

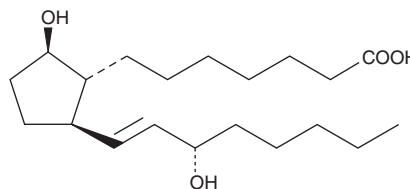
PRODUCT INFORMATION



11-deoxy Prostaglandin F_{1β}

Item No. 15510

CAS Registry No.: 37785-99-2
Formal Name: 9β,15S-dihydroxy-prost-13E-en-1-oic acid
Synonyms: 9β,11-deoxy PGF_{1α}, 11-deoxy PGF_{1β}
MF: C₂₀H₃₆O₄
FW: 340.5
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 F_{1β} (11-deoxy PGF_{1β}) is supplied as a crystalline solid. A stock solution may be made by dissolving the 11-deoxy PGF_{1β} in the solvent of choice, which should be purged with an inert gas. 11-deoxy PGF_{1β} is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 11-deoxy PGF_{1β} in ethanol is approximately 75 mg/ml and approximately 50 mg/ml in DMSO and DMF.

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 PGF_{1β} can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 11-deoxy PGF_{1β} 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 PGF_{1β} is a synthetic analog of PGF_{1β}. In contrast to PGF_{2α} and PGF_{1α}, 11-deoxy PGF_{1β} exhibits vasodepressor and bronchodilator activities in guinea pigs at a dose of 500 μg/kg.¹

Reference

1. Hall, D.W.R. and Jaitly, K.D. Structure-activity relationships in a series of 11-deoxy prostaglandins. *Prostaglandins* **11(3)**, 573-587 (1976).

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, 02/01/2024

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