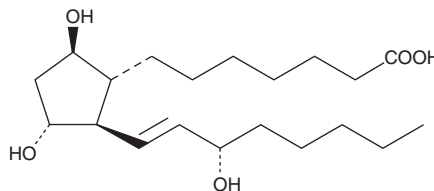


PRODUCT INFORMATION



Prostaglandin F_{1β} Item No. 15410

CAS Registry No.: 10164-73-5
Formal Name: 9β,11α,15S-trihydroxy-prost-13E-en-1-oic acid
Synonyms: 9β-PGF_{1α}, PGF_{1β}
MF: C₂₀H₃₆O₅
FW: 356.5
Purity: ≥98%
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

Prostaglandin F_{1β} (PGF_{1β}) is supplied as a crystalline solid. A stock solution may be made by dissolving the PGF_{1β} in the solvent of choice, which should be purged with an inert gas. PGF_{1β} is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 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 PGF_{1β} can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 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

PGF_{1β} is the C-9 epimer of PGF_{1α}. It was shown to enhance respiratory rate in experimental animals when administered intravenously.¹

Reference

1. Brookes, L.G. and Marshall, R.C. The effects of some prostaglandins on respiration in the rabbit. *J. Pharm. Pharmacol.* **26(Suppl)**, 80P-81P (1974).

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