

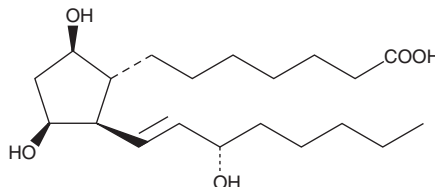
PRODUCT INFORMATION



11 β -Prostaglandin F_{1 β}

Item No. 15450

CAS Registry No.: 37785-86-7
Formal Name: 9 β ,11 β ,15S-trihydroxy-prost-13E-en-1-oic acid
Synonyms: 9 β ,11 β -PGF_{1 α} , 11 β -PGF_{1 β} , 11-*epi* PGF_{1 β}
MF: C₂₀H₃₆O₅
FW: 356.5
Purity: \geq 99%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

11 β -Prostaglandin F_{1 β} is supplied as a crystalline solid. A stock solution may be made by dissolving the 11 β -prostaglandin F_{1 β} in the solvent of choice. 11 β -Prostaglandin F_{1 β} is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 11 β -prostaglandin F_{1 β} 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 β -prostaglandin F_{1 β} can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 11 β -prostaglandin F_{1 β} 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 β -PGF_{1 β} is the stereoisomer of PGF_{1 α} with both C-9 and C-11 hydroxyls inverted.

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, 11/20/2018

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