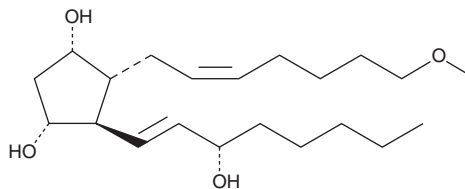


PRODUCT INFORMATION



Prostaglandin F_{2α} Alcohol methyl ether Item No. 16014

CAS Registry No.: 143656-18-2
Formal Name: 1-methoxy-prosta-5Z,13E-diene-9α,11α,15S-triol
Synonyms: AGN 191129, PGF_{2α}-OMe
MF: C₂₁H₃₈O₄
FW: 354.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Prostaglandin F_{2α} alcohol methyl ether (PGF_{2α}-OMe) 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 PGF_{2α}-OMe in these solvents is approximately 9, 15, and 7 mg/ml, respectively.

PGF_{2α}-OMe is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of PGF_{2α}-OMe should be diluted with the aqueous buffer of choice. The solubility of PGF_{2α}-OMe in PBS (pH 7.2) is approximately 0.16 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGF_{2α}-OMe is an analog of PGF_{2α} in which the C-1 carboxyl group has been replaced an O-methyl ether. The compound is reported to retain ocular hypotensive properties, but the receptors which mediate this activity have not been clearly documented.^{1,2}

References

1. Woodward, D.F. and Chan, M.F. Preparation of PGF-1 alcohols as ocular hypotensives. *Allergen, Inc.* **5,238,961** (1993).
2. Woodward, D.F., Krauss, A.H.P., Chen, J., *et al.* Replacement of the carboxylic acid group of prostaglandin F_{2α} with a hydroxyl or methoxy substituent provides biologically unique compounds. *Br. J. Pharmacol.* **130(8)**, 1933-1943 (2000).

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, 08/11/2022

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