

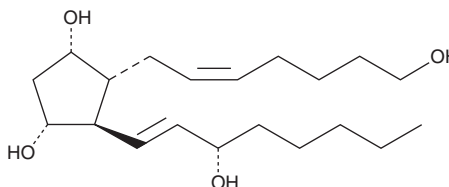
PRODUCT INFORMATION



Prostaglandin F_{2α} Alcohol

Item No. 16012

CAS Registry No.: 13261-27-3
Formal Name: prosta-5Z,13E-diene-1,9α,11α,15S-tetrol
Synonyms: AGN 190190, PGF_{2α}-OH, ZK 62717
MF: C₂₀H₃₆O₄
FW: 340.5
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 (PGF_{2α}-OH) 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 (DMF) purged with an inert gas can be used. The solubility of PGF_{2α}-OH in ethanol is approximately 9 mg/ml and approximately 29 mg/ml in DMSO and DMF.

PGF_{2α}-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of PGF_{2α}-OH should be diluted with the aqueous buffer of choice. The solubility of PGF_{2α}-OH in PBS (pH 7.2) is approximately 0.36 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PGF_{2α}-OH is an analog of PGF_{2α} in which the C-1 carboxyl group has been reduced to a primary alcohol. Biological studies on rat uterus show it to be a weak agonist.¹ The compound is reported to retain ocular hypotensive properties,² but the nature of the receptors which mediate these effects is disputed.³

References

1. Brambaifa, N. Luteolytic potency of PGD and PGF_{2α} derivatives. *Eur. J. Obstet. Gynecol. Reprod. Biol.* **27(4)**, 319-325 (1988).
2. Woodward, D.F. and Chan, M.F. Preparation of PGF-1 alcohols as ocular hypotensives. *Allergen, Inc.* **5,238,961** (1993).
3. 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, 08/10/2022

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