

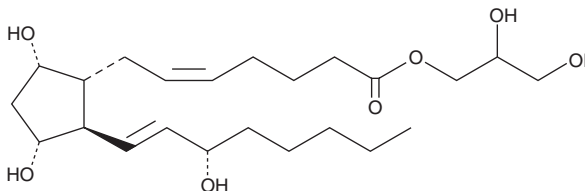
PRODUCT INFORMATION



Prostaglandin F_{2α}-1-glyceryl ester

Item No. 10139

CAS Registry No.: 43042-79-1
Formal Name: 9α,11α,15S-trihydroxy-prosta-5Z,13E-dien-1-oic acid, -1-glyceryl ester
Synonym: PGF_{2α}-1-glyceryl ester
MF: C₂₃H₄₀O₇
FW: 428.6
Purity: ≥98% (as a 9:1 mixture of the 1- and 2- glyceryl esters)
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that prostaglandin F_{2α}-1-glyceryl ester (PGF_{2α}-1-glyceryl ester) be stored as supplied at -20°C. It should be stable for at least one year.

PGF_{2α}-1-glyceryl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of PGF_{2α}-1-glyceryl ester in these solvents is approximately 20 mg/ml.

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. If an organic solvent-free solution of PGF_{2α}-1-glyceryl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of PGF_{2α}-1-glyceryl ester in PBS (pH 7.2) is approximately 500 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-Arachidonoyl glycerol (2-AG; Item No. 62160) has been isolated from porcine brain,¹ and has been characterized as the natural endocannabinoid ligand for the central cannabinoid receptor.² Incubation of 2-AG with cyclooxygenase-2 and specific PGH₂ isomerases *in vitro* in cell culture and isolated enzyme preparations results in PG glyceryl ester formation.³ The biosynthesis of PGH, PGD, PGE, PGF, and thromboxane A-2-glyceryl ester compounds have all been documented. The 2-glyceryl ester moiety equilibrates rapidly (within minutes) with the more stable 1-glyceryl ester, producing a 10:90 2:-1-glyceryl ester mixture in typical aqueous media. While the stability and metabolism of PGF_{2α}-1-glyceryl ester has been investigated, little is known about its intrinsic biological activity.⁴

References

1. Sugiura, T., Kodaka, T., Kondo, S., *et al. Biochem. Biophys. Res. Commun.* **229**, 58-64 (1996).
2. Sugiura, T., Kodaka, T., Kondo, S., *et al. J. Biochem.* **122**, 890-895 (1997).
3. Kozak, K.R., Crews, B.C., Morrow, J.D., *et al. J. Biol. Chem.* **277(47)**, 44877-44885 (2002).
4. Kozak, K.R., Crews, B.C., Ray, J.L., *et al. J. Biol. Chem.* **276(40)**, 36993-36998 (2001).

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, 08/13/2015

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