

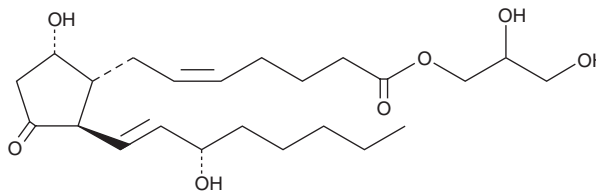
PRODUCT INFORMATION



Prostaglandin D₂-1-glyceryl ester

Item No. 12015

CAS Registry No.: 309260-52-4
Formal Name: 9 α ,15S-dihydroxy-11-oxo-prosta-5Z,13E-dien-1-oic acid, 1-glyceryl ester
Synonym: PGD₂-1-glyceryl ester
MF: C₂₃H₃₈O₇
FW: 426.6
Purity: \geq 95% (as a 9:1 mixture of the 1- and 2-glyceryl esters)
Stability: \geq 1 year at -80°C
Supplied as: A solution in acetonitrile



Laboratory Procedures

For long term storage, we suggest that prostaglandin D₂-1-glyceryl ester (PGD₂-1-glyceryl ester) be stored as supplied at -80°C. It should be stable for at least one year.

PGD₂-1-glyceryl ester is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile 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 PGD₂-1-glyceryl ester in these solvents is approximately 10 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 PGD₂-1-glyceryl ester is needed, it can be prepared by evaporating the acetonitrile and directly dissolving the neat oil in aqueous buffers. The solubility of PGD₂-1-glyceryl ester in PBS (pH 7.2) is approximately 0.1 mg/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, and has been characterized as the natural endocannabinoid ligand for the CB₁ receptor.^{1,2} Incubation of 2-AG with COX-2 and specific prostaglandin H₂ (PGH₂; Item No. 17020) isomerases in cell cultures and isolated enzyme preparations results in prostaglandin glycerol ester formation.³ The biosynthesis of PGH, PGD, PGE, PGF, and TXA-2-glyceryl ester compounds have all been documented. In RAW 264.7 cells, PGD₂-2-glyceryl ester is the main COX metabolite.³ The 2-glyceryl ester moiety equilibrates rapidly (within minutes) with the more stable 1-glyceryl ester, producing a 10:90 mixture of the 1- and 2-glyceryl esters in typical aqueous media. While the stability and metabolism of these PG products have been investigated, little is known about their 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, 03/03/2016

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