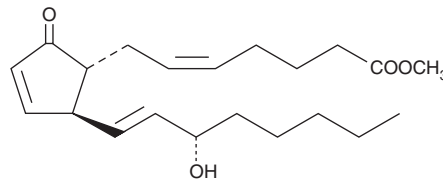


PRODUCT INFORMATION



Prostaglandin A₂ methyl ester Item No. 9000185

CAS Registry No.: 31753-19-2
Formal Name: 15S-hydroxy-9-oxo-prosta-5Z,10,13E-trien-1-oic acid, methyl ester
Synonyms: Medullin methyl ester, PGA₂ methyl ester
MF: C₂₁H₃₂O₄
FW: 348.5
Purity: ≥98%
UV/Vis.: λ_{max}: 215 nm
Supplied as: A solution in ethanol
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 A₂ methyl ester (PGA₂ methyl 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 PGA₂ methyl ester in these solvents is approximately 50, and 75 mg/ml, respectively.

PGA₂ methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of PGA₂ methyl ester should be diluted with the aqueous buffer of choice. The solubility of PGA₂ methyl ester in PBS (pH 7.2) is approximately 2 mg/ml. Avoid adding PGA₂ to basic solutions (pH > 7.4), since base treatment will degrade the PGA₂ to PGB₂. We do not recommend storing the aqueous solution for more than one day.

Description

Prostaglandin A₂ (PGA₂) methyl ester is a methylated derivative of the prostaglandin PGA₂ (Item No. 10210).¹ PGA₂ methyl ester has been found in *P. homomalla*.

Reference

- Schneider, W.P., Bundy, G.L., Lincoln, F.H., *et al.* Isolation and chemical conversions of prostaglandins from *Plexaura homomalla*: Preparation of prostaglandin E₂, prostaglandin F_{2α}, and their 5,6-trans isomers. *J. Am. Chem. Soc.* **99**(4), 1222-1232 (1977).

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, 02/07/2024

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