

PRODUCT INFORMATION



15(S)-acetate Prostaglandin A₂ methyl ester

Item No. 28195

CAS Registry No.: 36323-03-2
Formal Name: (5Z,13E,15S)-15-(acetyloxy)-9-oxo-prosta-5,10,13-trien-1-oic acid, methyl ester

Synonyms: 15-acetyl PGA₂ methyl ester, 15S-O-acetyl-PGA₂ methyl ester, 15S-PGA₂ methyl ester 15-acetate

MF: C₂₃H₃₄O₅

FW: 390.5

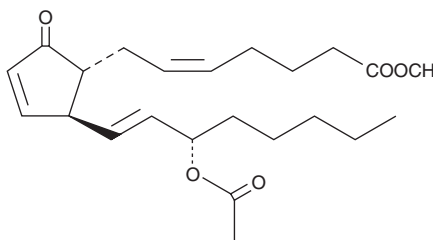
Purity: ≥95%

UV/Vis.: λ_{max}: 216 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

15(S)-acetate Prostaglandin A₂ 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 15(S)-acetate prostaglandin A₂ methyl ester in these solvents is approximately 30 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 15(S)-acetate prostaglandin A₂ methyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15(S)-acetate prostaglandin A₂ methyl ester in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15(S)-acetate Prostaglandin A₂ methyl ester is a 15-acetylated derivative of prostaglandin A₂ methyl ester (Item No. 9000185). It has been found in *P. homomalla*.¹

Reference

1. Ciereszko, L.S., Gopichand, Y., Schmitz, F.J., *et al.* Prostaglandin (15S)-PGA₂ derivatives in the gorgonian *Plexaura homomalla* (Esper), forma *kükenthali* Moser. *Experientia* **41(1)**, 37-38 (1985).

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, 01/31/2024

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