

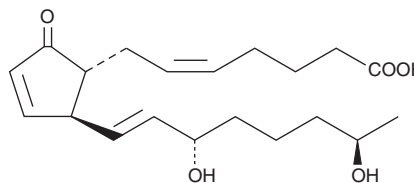
PRODUCT INFORMATION



19(R)-hydroxy Prostaglandin A₂

Item No. 10295

CAS Registry No.: 52087-58-8
Formal Name: 9-oxo-15S,19R-dihydroxy-prosta-5Z,10,13E-trien-1-oic acid
MF: C₂₀H₃₀O₅
FW: 350.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol
UV/Vis.: λ_{max}: 217 nm



Laboratory Procedures

For long term storage, we suggest that 19(R)-hydroxy Prostaglandin A₂ (19(R)-hydroxy PGA₂) be stored as supplied at -20°C. It should be stable for at least one year.

19(R)-hydroxy PGA₂ 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 19(R)-hydroxy PGA₂ in these solvents is approximately 50 and 75 mg/ml, respectively.

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 aqueous solution of 19(R)-hydroxy PGA₂ is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 19(R)-hydroxy PGA₂ in PBS (pH 7.2) is approximately 2.4 mg/ml. Store aqueous solutions of 19(R)-hydroxy PGA₂ on ice and use within 12 hours of preparation.

Description

19(R)-hydroxy PGA₂ has been found in human seminal plasma. Non-enzymatic dehydration of the 19(R)-hydroxy PGE₂, also found in semen, is presumed to be the source of this metabolite.¹

Reference

1. Hamberg, M. and Samuelsson, B. Prostaglandins in human seminal plasma. *J. Biol. Chem.* **241**, 257-263 (1966).

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, 08/18/2015

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