

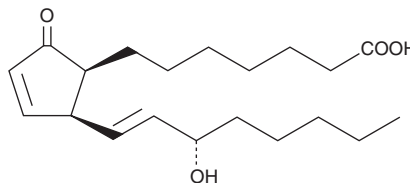
PRODUCT INFORMATION



8-iso Prostaglandin A₁

Item No. 10035

CAS Registry No.: 211186-29-7
Formal Name: 9-oxo-15S-hydroxy-(8β)-prosta-10,13E-dien-1-oic acid
Synonyms: 8-*epi* PGA₁, 8-*iso* PGA₁
MF: C₂₀H₃₂O₄
FW: 336.5
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A solution in methyl acetate
UV/Vis.: λ_{max}: 217 nm



Laboratory Procedures

For long term storage, we suggest that 8-*iso* prostaglandin A₁ (8-*iso* PGA₁) be stored as supplied at -20°C. It should be stable for at least two years.

8-*iso* PGA₁ is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate 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 8-*iso* PGA₁ in these solvents is approximately 100, 50, and 75 mg/ml, respectively. 8-*iso* PGA₁ will be stable for at least six months in these solvents if stored at -20°C.

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 8-*iso* PGA₁ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 8-*iso* PGA₁ in PBS, pH 7.2, is approximately 2.4 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

8-*iso* PGA₁ is an isoprostane and a member in a large family of prostanoids of non-cyclooxygenase origin. It occurs as a common minor impurity in most commercial preparations of PGE₁. The biological activity of 8-*iso* PGA₁ has not been studied in depth or reported in the literature.

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, 07/30/2015

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