

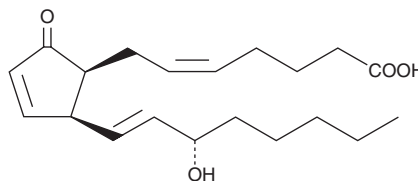
PRODUCT INFORMATION



8-iso Prostaglandin A₂

Item No. 10235

CAS Registry No.: 474391-66-7
Formal Name: 9-oxo-15S-hydroxy-(8β)-prosta-5Z,10,13E-trien-1-oic acid
Synonym: 8-*epi* PGA₂
MF: C₂₀H₃₀O₄
FW: 334.5
Purity: ≥98%
Stability: ≥1 year 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 A₂ (8-*iso* PGA₂) be stored as supplied at -20°C. It should be stable for at least one year.

8-*iso* PGA₂ 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* PGA₂ in these solvents is approximately 100, 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 solution of 8-*iso* PGA₂ 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* PGA₂ 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

Isoprostanes are PG-like compounds produced *in vivo* by free radical-catalyzed peroxidation of arachidonoyl-containing lipids. 8-*iso* PGA₂ is one of several isoprostanes produced from peroxidation of arachidonic acid esterified in phospholipids.¹ 8-*iso* PGA₂ is the dehydration product of 8-*iso* PGE₂, a potent renal vasoconstrictor.² Evidence for the *in vivo* production of 8-*iso* PGA₂ has been shown in rat liver under conditions of oxidative stress.³ There are no published studies of the pharmacological properties of 8-*iso* PGA₂.

References

1. Morrow, J.D., Minton, T.A., Mukundan, C.R., *et al.* Free radical-induced generation of isoprostanes *in vivo*. Evidence for the formation of D-ring and E-ring isoprostanes. *J. Biol. Chem.* **269**, 4317-4326 (1994).
2. Longmire, A.W., Roberts, L.J., and Morrow, J.D. Actions of the E₂-isoprostane, 8-*iso*-PGE₂, on the platelet thromboxane/endoperoxide receptor in humans and rats: Additional evidence for the existence of a unique isoprostane receptor. *Prostaglandins* **48**, 247-256 (1994).
3. Chen, Y., Zackert, W.E., Robers, L.J.II., *et al.* Evidence for the formation of a novel cyclopentenone isoprostane, 15-A_{2t}-isoprostane (8-*iso*-prostaglandin A₂) *in vivo*. *Biochem. Biophys. Acta* **1436**, 550-556 (1999).

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, 08/13/2015

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