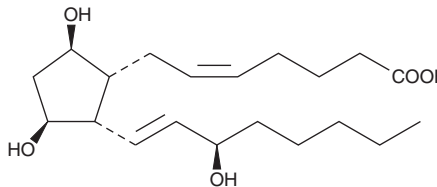


PRODUCT INFORMATION



ent-8-*iso* Prostaglandin F_{2a} Item No. 10011545

CAS Registry No.: 159812-83-6
Formal Name: 9*b*,11*b*,15*R*-trihydroxy-(12*a*)-prosta-5*Z*,13*E*-dien-1-oic acid
Synonyms: *ent*-8-*epi* PGF_{2a}, *ent*-8-*iso* PGF_{2a}, *ent*-15-F_{2t}-Isoprostane
MF: C₂₀H₃₄O₅
FW: 354.4
Purity: ≥98%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ent-8-*iso* PGF_{2a} is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile 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 *ent*-8-*iso* PGF_{2a} in these solvents is approximately 100 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 *ent*-8-*iso* PGF_{2a} is needed, it can be prepared by evaporating the acetonitrile and directly dissolving the neat oil in aqueous buffers. The solubility of *ent*-8-*iso* PGF_{2a} in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Isoprostanes are produced by the non-enzymatic, free radical peroxidation of arachidonic acid. They have been used as biomarkers of oxidative stress, but also have been found to have a potent biological activity. *ent*-8-*iso* PGF_{2a} is a potent vasoconstrictor of porcine retinal and brain microvessels with EC₅₀ values of 31 and 54 nM, respectively.¹

Reference

- Hou, X., Robers, L.J.II., Gobeil, F., Jr., *et al.* Isomer-specific contractile effects of a series of synthetic F₂-isoprostanes on retinal and cerebral microvasculature. *Free Radic. Biol. Med.* **36(2)**, 163-172 (2004).

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, 05/10/2024

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