

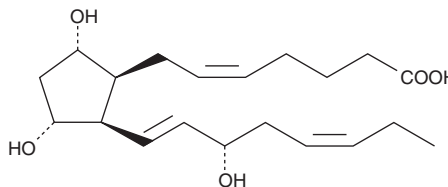
PRODUCT INFORMATION



8-iso Prostaglandin F_{3α}

Item No. 16992

CAS Registry No.: 7045-31-0
Formal Name: (8β)-9α,11α,15S-trihydroxy-prosta-5Z,13E,17Z-trien-1-oic acid
Synonyms: 8-*epi* PGF_{3α}, 8-*iso* PGF_{3α}
MF: C₂₀H₃₂O₅
FW: 352.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

8-*iso* Prostaglandin F_{3α} (8-*iso* PGF_{3α}) 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* PGF_{3α} in these solvents is approximately 100 mg/ml. 8-*iso* PGF_{3α} is soluble in 10 mM Na₂CO₃ at a concentration of approximately 6.5 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 8-*iso* PGF_{3α} 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* PGF_{3α} in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

8-*iso* PGF_{3α} is an isoprostane produced from the free-radical peroxidation of EPA. Little is known about the biological activity of 8-*iso* PGF_{3α}. There is one report that it is inactive in a TP receptor mediated assay of human platelet shape change, where 8-*iso* PGF_{2α} has an ED₅₀ value of 1 μM.¹

Reference

1. Praticò, D., Smith, E.M., Violi, F., *et al.* Local amplification of platelet function by 8-*epi* prostaglandin F_{2α} is not mediated by thromboxane receptor isoforms. *J. Biol. Chem.* **271**(25), 14916-14924 (1996).

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, 03/18/2024

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