

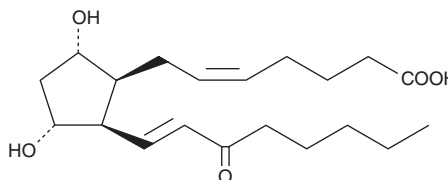
PRODUCT INFORMATION



8-iso-15-keto Prostaglandin F_{2α}

Item No. 16390

CAS Registry No.: 191919-01-4
Formal Name: 9α,11α-dihydroxy-15-oxo-(8β)-prosta-5Z,13E-dien-1-oic acid
Synonyms: 8-*epi*-15-keto PGF_{2α}, 8-*iso*-15-keto PGF_{2α}
MF: C₂₀H₃₂O₅
FW: 352.5
Purity: ≥95%
UV/Vis.: λ_{max}: 230 nm
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*-15-keto Prostaglandin F_{2α} (8-*iso*-15-keto PGF_{2α}) 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*-15-keto PGF_{2α} in these solvents is approximately 100 mg/ml. 8-*iso*-15-keto PGF_{2α} 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*-15-keto PGF_{2α} 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*-15-keto PGF_{2α} 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*-15-keto PGF_{2α} is a metabolite of the isoprostane 8-*iso* PGF_{2α} in rabbits, monkeys, and humans. 8-*iso* prostane (8-*iso* PGF_{2α}) is a prostaglandin-like product of non-specific lipid peroxidation.¹ In both humans and monkeys, exogenously infused 8-*iso* PGF_{2α} is converted primarily to metabolites having 2 or 4 carbon atoms removed from the top side chain by β-oxidation.² A similar pattern is observed when tritiated 8-*iso* PGF_{2α} is infused into rabbits.³ Early in the infusion (within 1-2 minutes) 8-*iso*-15-keto PGF_{2α} was a major component of the metabolite profile, which was comprised mostly of unmetabolized 8-*iso* PGF_{2α}. 8-*iso*-15-keto PGF_{2α} is a vasoconstrictor when tested on the rat isolated thoracic aorta, acting *via* the TP (thromboxane) receptor.⁴

Reference

1. Morrow, J.D., Hill, K.E., Burk, R.F., *et al.* A series of prostaglandin F₂-like compounds are produced *in vivo* in humans by a non-cyclooxygenase, free radical-catalyzed mechanism. *Proc. Natl. Acad. Sci. USA* **87(23)**, 9383-9387 (1990).
2. Chiabrando, C., Valagussa, A., Rivalta, C., *et al.* Identification and measurement of endogenous β-oxidation metabolites of 8-*epi*-prostaglandin F_{2α}. *J. Biol. Chem.* **274(3)**, 1313-1319 (1999).
3. Basu, S. Metabolism of 8-*iso*-prostaglandin F_{2α}. *FEBS Lett.* **428(112)**, 32-36 (1998).
4. Cracowski, J.L., Camus, L., Durand, T., *et al.* Response of rat thoracic aorta to F₂-isoprostane metabolites. *J. Cardiovasc. Pharmacol.* **39(3)**, 396-403 (2002).

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, 03/14/2024

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