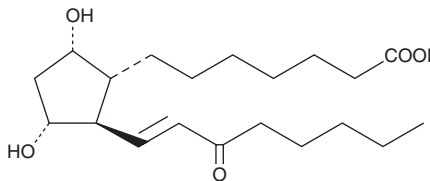


PRODUCT INFORMATION



15-keto Prostaglandin F_{1α} Item No. 15710

CAS Registry No.: 21562-58-3
Formal Name: 9α,11α-dihydroxy-15-oxo-prost-13E-en-1-oic acid
Synonym: 15-keto PGF_{1α}
MF: C₂₀H₃₄O₅
FW: 354.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232 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

15-keto Prostaglandin F_{1α} (15-keto PGF_{1α}) 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 (DMF) purged with an inert gas can be used. The solubility of 15-keto PGF_{1α} in ethanol is approximately 75 mg/ml and approximately 50 mg/ml in DMSO and DMF.

15-keto PGF_{1α} is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 15-keto PGF_{1α} should be diluted with the aqueous buffer of choice. The solubility of 15-keto PGF_{1α} in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15-keto PGF_{1α} is the initial metabolite of PGF_{1α} via 15-hydroxy PGDH. In mammals, oxidation of C-15 markedly attenuates receptor binding and activity. In fish, the 15-keto compounds serve as post-ovulatory pheromones and are more active than the parent prostaglandins.¹

Reference

1. Sorensen, P.W., Hara, T.J., Goetz, F.W., *et al.* F prostaglandins function as potent olfactory stimulants that comprise the postovulatory female sex pheromone in goldfish. *Biol. Reprod.* **39(5)**, 1039-1050 (1988).

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, 02/27/2023

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