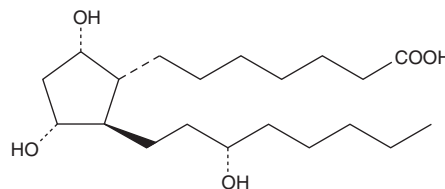


# PRODUCT INFORMATION



## 13,14-dihydro Prostaglandin F<sub>1α</sub> Item No. 15610

**CAS Registry No.:** 20592-20-5  
**Formal Name:** 9α,11α,15S-trihydroxy-prostan-1-oic acid  
**Synonyms:** PGF<sub>1α</sub>, 13,14-dihydro PGF<sub>1α</sub>  
**MF:** C<sub>20</sub>H<sub>38</sub>O<sub>5</sub>  
**FW:** 358.5  
**Purity:** ≥99%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

13,14-dihydro Prostaglandin F<sub>1α</sub> (13,14-dihydro PGF<sub>1α</sub>) is supplied as a crystalline solid. A stock solution may be made by dissolving the 13,14-dihydro PGF<sub>1α</sub> in the solvent of choice, which should be purged with an inert gas. 13,14-dihydro PGF<sub>1α</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 13,14-dihydro PGF<sub>1α</sub> in ethanol is approximately 75 mg/ml and approximately 50 mg/ml in DMSO and DMF.

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. Organic solvent-free aqueous solutions of 13,14-dihydro PGF<sub>1α</sub> can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 13,14-dihydro PGF<sub>1α</sub> in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

13,14-dihydro PGF<sub>1α</sub> is a potential metabolite of PGF<sub>1α</sub>.<sup>1</sup> Its biological activity has not been reported. The corresponding compound in the E<sub>1</sub> series (13,14-dihydro PGE<sub>1</sub>) retains biological activity similar to PGE<sub>1</sub>.<sup>2</sup>

### References

1. Ritzi, E.M., Boto, W.O., and Stylos, W.A. Measurement of initial prostaglandin F metabolites in medium of BALB/c 3T3 and SV3T3 mouse fibroblast cultures. *Biochem. Biophys. Res. Commun.* **63**(1), 179-186 (1975).
2. Ånggård, E. The biological activities of three metabolites of prostaglandin E<sub>1</sub>. *Acta Physiol. Scand.* **66**(4), 509-510 (1966).

#### 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, 02/06/2024

#### 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