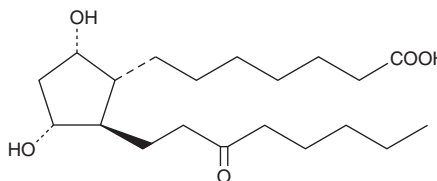


# PRODUCT INFORMATION



## 13,14-dihydro-15-keto Prostaglandin F<sub>1α</sub> Item No. 15670

**CAS Registry No.:** 29044-75-5  
**Formal Name:** 9α,11α-dihydroxy-15-oxo-prostan-1-oic acid  
**Synonym:** 13,14-dihydro-15-keto PGF<sub>1α</sub>  
**MF:** C<sub>20</sub>H<sub>36</sub>O<sub>5</sub>  
**FW:** 356.5  
**Purity:** ≥95%  
**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

13,14-dihydro-15-keto Prostaglandin F<sub>1α</sub> (13,14-dihydro-15-keto PGF<sub>1α</sub>) 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 13,14-dihydro-15-keto PGF<sub>1α</sub> in these solvents is approximately 50 mg/ml. 13,14-dihydro-15-keto PGF<sub>1α</sub> is stable for at least six months in these solvents if stored at -20°C.

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 13,14-dihydro-15-keto PGF<sub>1α</sub> is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 13,14-dihydro-15-keto 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-15-keto PGF<sub>1α</sub> is a metabolite of PGF<sub>1α</sub> that has been reported in the rat stomach.<sup>1</sup> The measurement of 13,14-dihydro-15-keto PGF<sub>1α</sub> can be used as a marker of the *in vivo* production of PGF<sub>1α</sub>.<sup>2</sup>

### References

1. Pace-Asciak, C.R., Morawska, K., and Wolfe, L.S. Metabolism of prostaglandin F<sub>1α</sub> by the rat stomach. *Biochim. Biophys. Acta* **218**, 288-295 (1970).
2. Sors, H., Maclouf, J., Pradelles, P., *et al.* The use of iodinated tracers for a sensitive radioimmunoassay of 13,14-dihydro-15-keto prostaglandin F<sub>α</sub>. *Biochim. Biophys. Acta* **486**, 553-564 (1977).

#### 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, 12/21/2018

#### 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