

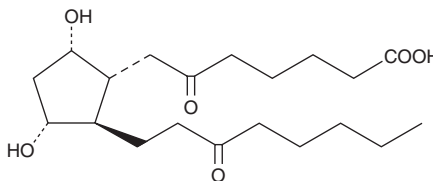
# PRODUCT INFORMATION



## 6,15-diketo-13,14-dihydro Prostaglandin F<sub>1α</sub>

Item No. 15270

**CAS Registry No.:** 63983-53-9  
**Formal Name:** 6,15-dioxo-9α,11α-dihydroxy-prostan-1-oic acid  
**Synonym:** 6,15-diketo-13,14-dihydro PGF<sub>1α</sub>  
**MF:** C<sub>20</sub>H<sub>34</sub>O<sub>6</sub>  
**FW:** 370.5  
**Purity:** ≥98% (mixture of tautomers)  
**Supplied as:** A 1 mg/ml 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

6,15-diketo-13,14-dihydro 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, or dimethyl formamide purged with an inert gas can be used. The solubility of 6,15-diketo-13,14-dihydro PGF<sub>1α</sub> in these solvents is approximately 16, 50, and 100 mg/ml, respectively.

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 aqueous solution of 6,15-diketo-13,14-dihydro 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 6,15-diketo-13,14-dihydro PGF<sub>1α</sub> in PBS (pH 7.2) is approximately 3.3 mg/ml. Store aqueous solutions of 6,15-diketo-13,14-dihydro PGF<sub>1α</sub> on ice and use within 12 hours of preparation.

### Description

6,15-diketo-13,14-dihydro PGF<sub>1α</sub> is a metabolite of PGI<sub>2</sub>. It was shown to enhance intracellular cAMP and cholesterol catabolism in bovine arterial smooth muscle cells.<sup>1</sup>

### Reference

1. Etingin, O.R., Weksler, B.B., and Hajjar, D.P. Cholesterol metabolism is altered by hydrolytic metabolites of prostacyclin in arterial smooth muscle cells. *J. Lipid Res.* **27**, 530-536 (1986).

#### 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, 12/16/2025

#### 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](http://WWW.CAYMANCHEM.COM)