

# PRODUCT INFORMATION



## 9-keto Fluprostenol

Item No. 16781

**CAS Registry No.:** 156406-33-6  
**Formal Name:** (5Z)-7-[(1R,2R,3R)-3-hydroxy-2-[(1E,3R)-3-hydroxy-4-[3-(trifluoromethyl)phenoxy]-1-buten-1-yl]-5-oxocyclopentyl]-5-heptenoic acid

**Synonym:** Fluprostenol Prostaglandin E<sub>2</sub>

**MF:** C<sub>23</sub>H<sub>27</sub>F<sub>3</sub>O<sub>6</sub>

**FW:** 456.5

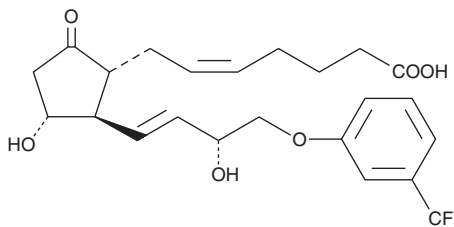
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 222, 277 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

9-keto Fluprostenol 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. 9-keto Fluprostenol is miscible in ethanol.

9-keto Fluprostenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 9-keto fluprostenol should be diluted with the aqueous buffer of choice. The solubility of 9-keto fluprostenol in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

9-keto Fluprostenol is an analog of prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) with structural modifications intended to give it a prolonged half-life and greater potency. Fluprostenol is a well-studied, potent analog of PGF<sub>2α</sub> and acts primarily through the FP receptor.<sup>1</sup> Oxidation at C-9 of fluprostenol yields 9-keto fluprostenol. It is anticipated that this analog will have strong affinity for EP receptors and act as a PGE<sub>2</sub> agonist.

### Reference

1. Dukes, M., Russell, W., and Walpole, A.L. Potent luteolytic agents related to prostaglandin F<sub>2α</sub>. *Nature* **250(464)**, 330-331 (1974).

#### 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, 12/01/2023

#### 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