

PRODUCT INFORMATION



(+)-Cloprostenoil isopropyl ester

Item No. 10010016

CAS Registry No.: 157283-66-4

Formal Name: (+)-9 α ,11 α ,15R-trihydroxy-16-(3-chlorophenoxy)-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-oic acid, isopropyl ester

Synonyms: (+)-5-cis Cloprostenoil isopropyl ester, (+)-16-m-chlorophenoxy tetranor Prostaglandin F_{2 α} isopropyl ester

MF: C₂₅H₃₅ClO₆

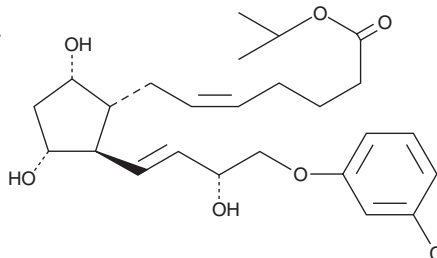
FW: 467.0

Purity: \geq 98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: \geq 2 years



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

Laboratory Procedures

(+)-Cloprostenoil isopropyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of (+)-cloprostenoil isopropyl ester in these solvents is approximately 100 mg/ml.

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 (+)-cloprostenoil isopropyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (+)-cloprostenoil isopropyl ester in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Cloprostenoil is a synthetic analog of prostaglandin F_{2 α} (PGF_{2 α}). It is an FP receptor agonist and a potent luteolytic agent in rats and hamsters. (+)-Cloprostenoil is the optically active, 15(R) enantiomer of cloprostenoil responsible for the majority of its biological activity. It is 200 times more potent than PGF_{2 α} in terminating pregnancy when given subcutaneously at a daily dose of 0.125 μ g/kg in rats and hamsters, without the side effects associated with PGF_{2 α} .¹ (+)-Cloprostenoil was also shown to be a potent inhibitor of rat adipose precursor differentiation in primary cultures with an IC₅₀ value of 3 x 10⁻¹² M.² (+)-Cloprostenoil isopropyl ester is a more lipid soluble form of cloprostenoil.

References

1. Dukes, M., Russell, W., and Walpole, A.L. Potent luteolytic agents related to prostaglandin F_{2 α} . *Nature* **250(464)**, 330-331 (1974).
2. Serrero, G. and Lepak, N.M. Prostaglandin F_{2 α} receptor (FP receptor) agonists are potent adipose differentiation inhibitors for primary culture of adipocyte precursors in defined medium. *Biochem. Biophys. Res. Commun.* **233(1)**, 200-202 (1997).

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. It is the responsibility of the purchaser to determine suitability for other applications.

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, 04/21/2021

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