

PRODUCT INFORMATION



15-keto Cloprostenol isopropyl ester

Item No. 24929

Formal Name: isopropyl (Z)-7-((1R,2R,3R,5S)-2-((E)-4-(3-chlorophenoxy)-3-oxobut-1-en-1-yl)-3,5-dihydroxycyclopentyl)hept-5-enoate

MF: C₂₅H₃₃ClO₆

FW: 465.0

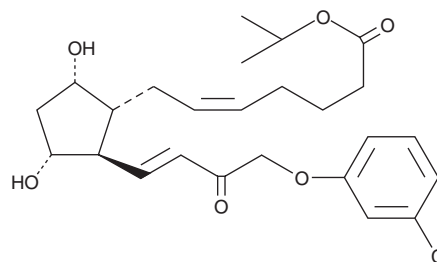
Purity: ≥98%

UV/Vis.: λ_{max}: 226 nm

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

15-keto Cloprostenol 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 15-keto cloprostenol 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 15-keto cloprostenol isopropyl ester is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15-keto cloprostenol 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

15-keto Cloprostenol isopropyl ester is a potential impurity found in commercial cloprostenol isopropyl ester preparations. (+)-Cloprostenol isopropyl ester (Item No. 10010016) is a more lipid soluble form of cloprostenol (Item No. 16764), a synthetic analog of prostaglandin F_{2α} (PGF_{2α}; Item No. 16010).

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, 01/31/2024

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