

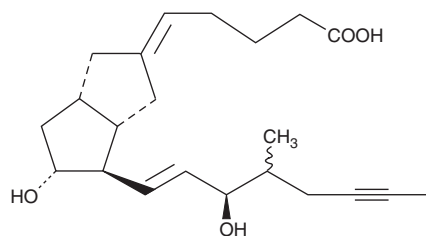
PRODUCT INFORMATION



5-*cis*-15(R)-Iloprost

Item No. 10008588

Formal Name:	6,9 α -methylene-11 α ,15R-dihydroxy-16-methyl-prosta-5Z,13E-dien-18-yn-1-oic acid
Synonyms:	5- <i>cis</i> -15(R)-Ciloprost
MF:	C ₂₂ H ₃₂ O ₄
FW:	360.5
Purity:	≥97%
Supplied as:	A 5 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

5-*cis*-15(R)-Iloprost 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 (DMF) purged with an inert gas can be used. The solubility of 5-*cis*-15(R)-iloprost in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

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 5-*cis*-15(R)-iloprost is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 5-*cis*-15(R)-iloprost in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-*cis*-15(R)-Iloprost is the C-5 *cis*-isomer and 15(R)-epimer of iloprost, a second generation structural analog of prostaglandin I₂ (prostacyclin). Iloprost displays ten-fold greater potency than the first generation stable prostacyclin analogs, typified by carbaprostacyclin.¹ There are no published studies of the pharmacological properties of 5-*cis*-15(R)-iloprost.

Reference

1. Schrör, K., Darius, H., Matzky, R., *et al.* The antiplatelet and cardiovascular actions of a new carbacyclin derivative (ZK 36 374) - equipotent to PGI₂ in vitro. *Naunyn Schmiedebergs Arch. Pharmacol.* **316(3)**, 252-255 (1981).

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, 01/15/2026

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