

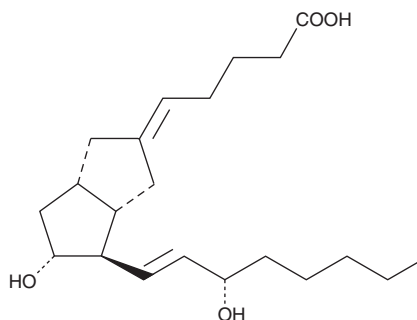
PRODUCT INFORMATION



5-cis Carbaprostacyclin

Item No. 18211

CAS Registry No.: 69609-77-4
Formal Name: (5Z)-5-[(3aS,4R,5R,6aS)-hexahydro-5-hydroxy-4-[(1E,3S)-3-hydroxy-1-octenyl]-2(1H)-pentalenyldiene]-pentanoic acid
Synonym: (5Z)-6a-Carbaprostaglandin I₂
MF: C₂₁H₃₄O₄
FW: 350.5
Purity: ≥99%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

5-cis Carbaprostacyclin is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-cis carbaprostacyclin in the solvent of choice, which should be purged with an inert gas. 5-cis Carbaprostacyclin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-cis carbaprostacyclin in these solvents is approximately 20, 5, and 10 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. Organic solvent-free aqueous solutions of 5-cis carbaprostacyclin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 5-cis carbaprostacyclin in PBS (pH 7.2) is approximately 0.08 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-cis Carbaprostacyclin is a stable analog of PGI₂ and an isomer of carbaprostacyclin. It is a weak inhibitor of human platelet aggregation with an IC₅₀ of 2.8 μM compared to 0.3 μM for carbaprostacyclin.^{1,2} 5-cis Carbaprostacyclin is a much weaker effector of rabbit mesenteric artery relaxation with an EC₅₀ of 104 μM compared to 5.9 μM for carbaprostacyclin. It even antagonizes the adenylate cyclase activation induced by carbaprostacyclin.²

References

1. Whittle, B.J.R., Moncada, S., Whiting, F., *et al.* Carbacyclin – a potent stable prostacyclin analogue for the inhibition of platelet aggregation. *Prostaglandins* **19(4)**, 605-627 (1980).
2. Corsini, A., Folco, G.C., Fumagalli, R., *et al.* (5Z)-carbacyclin discriminates between prostacyclin-receptors coupled to adenylate cyclase in vascular smooth muscle and platelets. *Br. J. Pharmacol.* **90(1)**, 255-261 (1987).

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, 02/15/2024

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