

PRODUCT INFORMATION



Iloprost-d₄

Item No. 9000366

Formal Name: 5-((3aS,4R,5R,6aS,E)-5-hydroxy-4-((3S,E)-3-hydroxy-4-methyloct-1-en-6-yn-1-yl)hexahydropentalen-2(1H)-ylidene)pentanoic-3,3,4,4-d₄ acid

Synonym: Ciloprost-d₄

MF: C₂₂H₂₈D₄O₄

FW: 364.5

Chemical Purity: ≥95% (Iloprost)

Deuterium

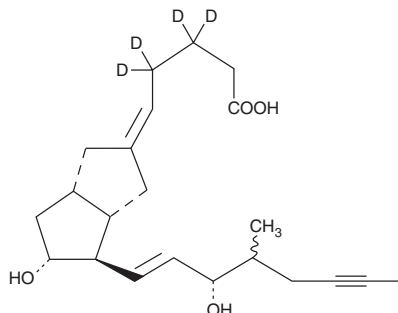
Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: ≥1 year

Special Conditions: Vasodilator



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

Laboratory Procedures

Iloprost-d₄ is intended for use as an internal standard for the quantification of iloprost (Item No. 18215) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Iloprost-d₄ 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 chloroform and dichloromethane purged with an inert gas can be used. Iloprost-d₄ is slightly soluble (0.1-1 mg/ml) in chloroform and dichloromethane.

Description

Iloprost is an agonist of the IP receptor and the prostaglandin E₂ (PGE₂) receptor subtype EP₁ and a derivative of PGI₂.^{1,2} It selectively binds to IP and EP₁ receptors (K_i = 11 nM for both) over EP₂, EP₄, DP, FP, and TP receptors (K_s = 1,870, 284, 1,035, 619, and 6,487 nM, respectively) but also binds to the EP₃ receptor (K_i = 56 nM).¹ Iloprost increases cAMP levels in HEK293 cells expressing IP or EP₃ receptors (EC₅₀s = 0.37 and 27.5 nM, respectively) and increases calcium levels in HEK293 cells expressing the EP₁ receptor (EC₅₀ = 0.3 nM).² It inhibits ADP-, thrombin-, and collagen-induced platelet aggregation in isolated human platelet-rich plasma (IC₅₀s = 1.07, 0.71, and 0.24 nM, respectively).³ Iloprost (100 ng/kg per minute) increases the time to occlusive coronary artery thrombosis in a porcine model of electrically induced coronary artery thrombosis.⁴ Aerosolized administration of iloprost (130-1,300 ng/kg per minute) reduces right ventricular systolic pressure and reverses vascular remodeling in rats in a model of chronic pulmonary hypertension induced by the alkaloid monocrotaline (Item No. 16666).⁵ Formulations containing iloprost have been used in the treatment of pulmonary arterial hypertension and severe frostbite.

References

1. Abramovitz, M., Adam, M., Boie, Y., et al. *Biochim. Biophys. Acta* **1483**(2), 285-293 (2000).
2. Whittle, B.J., Silverstein, A.M., Mottola, D.M., et al. *Biochem. Pharmacol.* **84**(1), 68-75 (2012).
3. Schrör, K., Darius, H., Matzky, R., et al. *Naunyn Schmiedeberg's Arch. Pharmacol.* **316**(3), 252-255 (1981).
4. van der Giessen, W.J., Mooi, W.J., Rutteman, A.M., et al. *Thromb. Res.* **36**(1), 45-51 (1984).
5. Schermuly, R.T., Yilmaz, H., Ghofrani, H.A., et al. *Am. J. Respir. Crit. Care Med.* **172**(3), 358-363 (2005).

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, 07/03/2024

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