

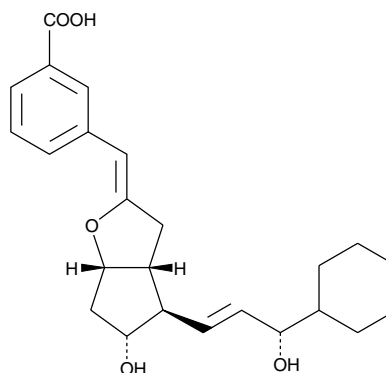
Product Information



Taprostene free acid

Item No. 10011348

CAS Registry No.: 108945-35-3
Formal Name: 3-[(Z)-[(3aR,4R,5R,6aS)-4-[(1E,3S)-3-cyclohexyl-3-hydroxy-1-propenyl]hexahydro-5-hydroxy-2H-cyclopenta[b]furan-2-ylidene)methyl]-benzoic acid
Synonyms: CG 4203, Rheocyclan
MF: C₂₄H₃₀O₅
FW: 398.5
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 233, 270 nm



Laboratory Procedures

For long term storage, we suggest that taprostene free acid be stored as supplied at -20°C. It should be stable for at least one year.

Taprostene free acid is supplied as a crystalline solid. For biological experiments, we suggest that organic solvent-free aqueous solutions of taprostene free acid be prepared by directly dissolving the crystalline compound in water. We do not recommend storing the aqueous solution for more than one day.

Taprostene is a stable prostacyclin (PGI₂) analog and agonist of the prostacyclin receptor, IP. It does not activate the PGE₂ receptor EP₄,¹ which, like IP, promotes vascular smooth muscle relaxation when stimulated. Taprostene has been used extensively to study the role of the IP receptor in tissue preparations and *in vivo*.¹⁻⁴ It has also been used in the screening and evaluation of potential IP antagonists.⁵

References

1. Jones, R.L. and Chan, K. Distinction between relaxations induced *via* prostanoid EP₄ and IP₁ receptors in pig and rabbit blood vessels. *Br. J. Pharmacol.* **134**, 313-324 (2001).
2. Tam, F.S.F., Chan, K., Bourreau, J.-P., *et al.* The mechanisms of enhancement and inhibition of field stimulation responses of guinea-pig vas deferens by prostacyclin analogues. *Br. J. Pharmacol.* **121**, 1413-1421 (1997).
3. Virgolini, I., Fitscha, P., Sinzinger, H., *et al.* Effects of taprostene, a stable prostacyclin analogue, on haemodynamics, platelet function and arachidonate metabolism in healthy volunteers. *Eur. J. Clin. Pharmacol.* **38**, 347-350 (1990).
4. Johnson, G., III, Furlan, L.E., Aoki, N., *et al.* Endothelium and myocardial protecting actions of taprostene, a stable prostacyclin analogue, after acute myocardial ischemia and reperfusion in cats. *Circ. Res.* **66**, 1362-1370 (1990).
5. Ayer, L.M., Wilson, S.M., Traves, S.L., *et al.* 4,5-dihydro-1H-imidazol-2-yl)-[4-(4-isopropoxy-benzyl)-phenyl]-amine (RO1138452) is a selective, pseudo-irreversible orthosteric antagonist at the prostacyclin (IP)-receptor expressed by human airway epithelial cells: IP-receptor-mediated inhibition of CXCL9 and CXCL10 release. *J. Pharmacol. Exp. Ther.* **324**(2), 815-826 (2008).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011348

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 06/24/2014

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com