

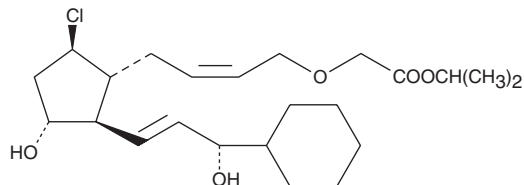
PRODUCT INFORMATION



ZK 118182 isopropyl ester

Item No. 13673

CAS Registry No.: 154927-31-8
Formal Name: [[[2Z]-4-[(1R,2R,3R,5R)-5-chloro-2-[(1E,3S)-3-cyclohexyl-3-hydroxy-1-propenyl]-3-hydroxycyclopentyl]-2-butenyl]oxy]-acetic acid, 1-methylethyl ester
MF: C₂₃H₃₇ClO₅
FW: 429.0
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

ZK 118182 isopropyl ester 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 purged with an inert gas can be used. The solubility of ZK 118182 isopropyl ester in these solvents is approximately 12.5, 2, and 3 mg/ml, respectively.

ZK 118182 isopropyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of ZK 118182 isopropyl ester should be diluted with the aqueous buffer of choice. ZK 118182 isopropyl ester has a solubility of 0.1 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Elevated intraocular pressure (IOP) is an important risk factor in developing glaucoma. Certain prostaglandins (PGs) such as PGF_{2α} and PGD₂, have been shown to reduce IOP. ZK 118182 is a PG analog that has potent DP-agonist activity (EC₅₀ = 16.5 nM) and a high nanomolar affinity for the DP receptor (K_i = 74 nM).¹ ZK 118182 isopropyl ester is a prodrug formulation of ZK 118182 designed to enhance corneal absorption. At 0.03 μg, ZK 118182 isopropyl ester has been shown to lower monkey and rabbit IOP 46% and 20%, respectively, two hours after dosing, demonstrating a much more potent effect compared to the free acid.²

References

1. Sharif, N.A., Williams, G.W., Crider, J.Y., *et al.* Molecular pharmacology of the DP/EP2 class prostaglandin AL-6598 and quantitative autoradiographic visualization of DP and EP2 receptor sites in human eyes. *J. Ocul. Pharmacol. Ther.* **20(6)**, 489-508 (2004).
2. Hellberg, M.R., Conrow, R.E., Sharif, N.A., *et al.* 3-Oxa-15-cyclohexyl prostaglandin DP receptor agonists as topical antiglaucoma agents. *Bioorg. Med. Chem.* **10(6)**, 2031-2049 (2002).

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, 03/08/2024

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