

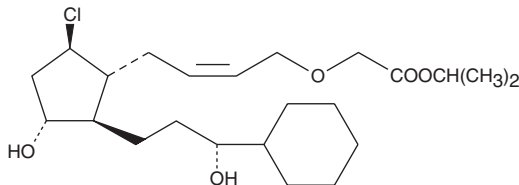
PRODUCT INFORMATION



AL 6598

Item No. 13620

CAS Registry No.: 170291-06-2
Formal Name: 2-[[[(2Z)-4-[(1R,2R,3R,5R)-5-chloro-2-[(3R)-3-cyclohexyl-3-hydroxypropyl]-3-hydroxycyclopentyl]-2-buten-1-yl]oxy]acetic acid, 1-methylethyl ester
MF: C₂₃H₃₉ClO₅
FW: 431.0
Purity: ≥97%
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

AL 6598 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 AL 6598 in ethanol is approximately 10 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

Elevated intraocular pressure (IOP) is an important risk factor in developing glaucoma. Certain prostaglandins such as PGF_{2α} and PGD₂, have been shown to reduce IOP. AL 6598 is the isopropyl ester prodrug of AL 6556, a PGD₂ receptor agonist that binds to DP receptors with a K_i value of 3.2 μM and demonstrates an EC₅₀ value of 0.80 μM in an *in vitro* functional assay.¹ Designed to enhance corneal absorption, AL 6598 produces a maximum 53% drop in IOP of the ocular hypertensive monkey with a 1 μg dose given twice daily.¹

Reference

1. 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, 04/13/2026

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