

PRODUCT INFORMATION



Latanoprost *tris*(triethylsilyl) ether

Item No. 37396

CAS Registry No.: 477884-78-9
Formal Name: (5Z)-7-[(1R,2R,3R,5S)-2-[(3R)-5-phenyl-3-[(triethylsilyl)oxy]pentyl]-3,5-bis[(triethylsilyl)oxy]cyclopentyl]-5-heptenoic acid, 1-methylethyl ester

Synonym: Latanoprost-related Compound D

MF: C₄₄H₈₂O₅Si₃

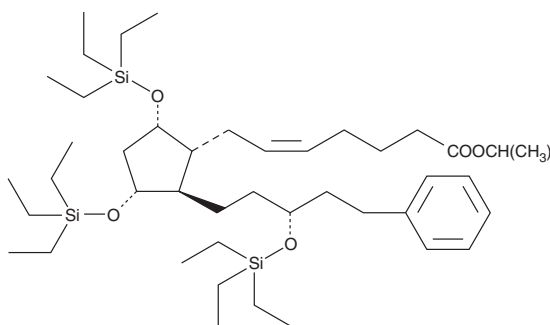
FW: 775.4

Purity: ≥95%

Supplied as: A neat oil

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Latanoprost *tris*(triethylsilyl) ether is supplied as a neat oil. A stock solution may be made by dissolving the latanoprost *tris*(triethylsilyl) ether in the solvent of choice, which should be purged with an inert gas. Latanoprost *tris*(triethylsilyl) ether is soluble in chloroform. Latanoprost *tris*(triethylsilyl) ether is slightly soluble in acetonitrile (warmed, sonicated).

Description

Latanoprost *tris*(triethylsilyl) ether is a precursor in the synthesis of the prostaglandin F_{2α} (PGF_{2α}) receptor (FP receptor) agonist latanoprost (Item No. 16812).¹

Reference

1. Greenwood, A.K., Mchattie, D., Thompson, D.G., *et al.* Process for the preparation of prostaglandins and analogues thereof. *Resolution Chemicals Limited*. **WO02/096898A2** (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, 11/01/2022

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