

PRODUCT INFORMATION



Latanoprost Analytical Reference Standard

Item No. 9000244

CAS Registry No.: 130209-82-4

Formal Name: 9 α ,11 α ,15R-trihydroxy-17-phenyl-18,19,20-trinor-prost-5Z-en-1-oic acid, isopropyl ester

Synonyms: 17-phenyl-13,14-dihydro trinor PGF_{2 α} isopropyl ester, 17-phenyl-13,14-dihydro trinor Prostaglandin F_{2 α} isopropyl ester

MF: C₂₆H₄₀O₅

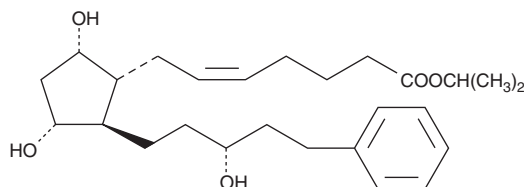
FW: 432.6

Purity: $\geq 99.0\%$

Supplied as: A neat oil

Storage: -20°C

Stability: ≥ 1 year



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

Description

Cayman's latanoprost analytical reference standard is a highly purified and characterized neat oil ideally suited as a reference material for assay determinations. Latanoprost is very hygroscopic. Limit atmospheric exposure and maintain the material in a tightly sealed container.

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, 01/21/2025E

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