

Latanoprost

Prostanoid receptor ligand

Prodrug (isopropyl ester) of 17-phenyl-13,14-dihydro-prostaglandin $F_{2\alpha}$.
The free acid form is a potent agonist at FP receptors. It is in clinical use as a topical antiglaucoma agent.

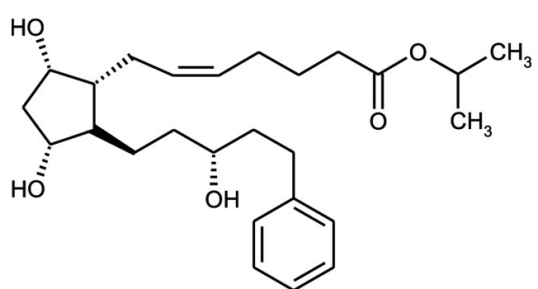
Ordering Information

[Order Online »](#)

BML-RA120-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Thick pale yellow oil.

CAS 130209-82-4

Couple Target Prostaglandin receptor, Prostanoid receptor

Couple Type Ligand

Formula $C_{26}H_{40}O_5$

Identity Determined by NMR.

MW 432.6

Purity ≥98% (Assay)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)