

Bimatoprost

FP receptor agonist

Stable analog of PGF₂α. Agonist at FP receptors (after *in vivo* hydrolysis to the free acid). It is an efficacious treatment for glaucoma. It also displays an unusual eye lash lengthening effect.

Citations: 2

[View Online »](#)

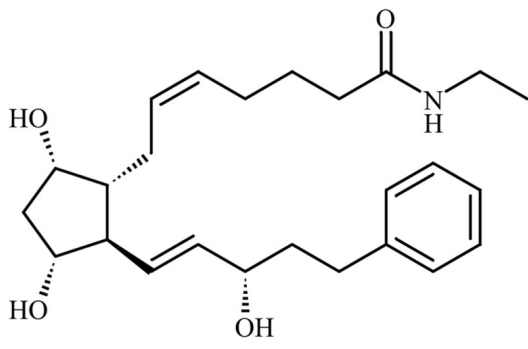
Ordering Information

[Order Online »](#)

BML-RA121-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(E,3S)-3-hydroxy-5-phenylpent-1-enyl]cyclopentyl]-N-ethylhept-5-enamide
Appearance	White to off-white powder.
CAS	155206-00-1
Couple Target	Prostanoid receptor
Couple Type	Activator, Ligand
Formula	$C_{25}H_{37}NO_4$
MW	415.6
Purity	≥97%
Solubility	Soluble in DMSO (25mg/ml) or ethanol (25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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