

D-erythro MAPP

Alkaline ceramidase inhibitor

A cell-permeable and specific inhibitor of alkaline ceramidase, which can raise intracellular ceramide levels by blocking the hydrolysis of ceramide to sphingosine and free fatty acid. Enhances the apoptotic response to ionizing radiation by enhancing the ceramide signal.

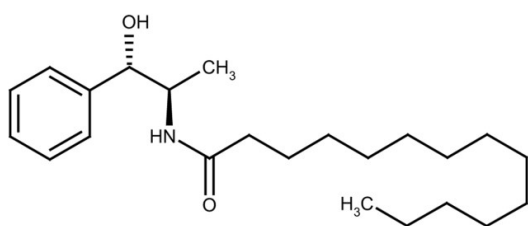
Ordering Information

[Order Online »](#)

BML-SL221-0005	5mg
BML-SL221-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(1S,2R-D-erythro-2-N-Myristoylamino)-1-phenyl-1-propanol
Appearance	White solid.
CAS	143492-38-0
Couple Target	Alkaline ceramidase
Couple Type	Inhibitor
Formula	$C_{23}H_{39}NO_2$
MW	361.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

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)