

N,N-Dimethylsphingosine, D-erythro

Sphingosine kinase inhibitor

Potent and specific inhibitor of sphingosine kinase ($IC_{50}=5\mu M$) which blocks conversion of sphingosine to sphingosine-1-phosphate (Prod. No. BML-SL140). Inhibitor of protein kinase C (PKC) which also stimulates Src-kinase. May affect expression of cell surface selectins, which in turn mediate leukocyte or tumor cell adhesion to endothelial cells and platelets. Induces apoptosis. Stereospecifically induces EGF receptor autophosphorylation producing EGF-like activity *in vitro* in the absence of EGF. Produced endogenously via ceramide catabolism. May be a target for neuropathic pain therapy.

Citations: 12

[View Online »](#)

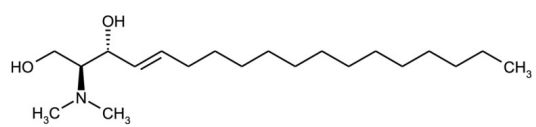
Ordering Information

[Order Online »](#)

BML-SL105-0025	25mg
----------------	------

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 at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N,N-Dimethyl-D-erythro-sphingosine
Appearance	White to off-white waxy solid or viscous oil.
CAS	119567-63-4
Couple Target	PKC, Sphingosine kinase
Couple Type	Inhibitor
Formula	$C_{20}H_{41}NO_2$
MW	327.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).
Source	Synthetic.
Technical Info / Product Notes	Note: Product is not sterile



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