

C6 Ceramide

Cell-permeable ceramide analog

A biologically active, cell-permeable, but nonphysiologic ceramide analog. It activates protein phosphatase 2A and MAP kinase (MAPK/ERK), suppresses insulin-induced tyrosine phosphorylation and inhibits glycoprotein traffic by the secretory pathway. It inhibits diacylglycerol accumulation and phospholipase D activation in fibroblasts. It induces an arrest at the G₀/G₁ transition of the cell cycle (10μM) and causes apoptosis (15μM or greater) in Molt-4 leukemia cells.

Citations: 9

[View Online »](#)

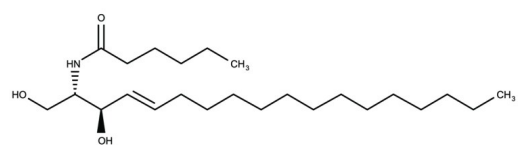
Ordering Information

[Order Online »](#)

BML-SL110-0005	5mg
BML-SL110-0025	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, as supplied, 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-Hexanoyl-D-erythro-sphingosine
Appearance	White waxy solid.
CAS	124753-97-5
Couple Target	Erk, MAPK, Serine/threonine-protein phosphatase
Couple Type	Activator
Formula	$C_{24}H_{47}NO_3$
MW	397.7
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (>25mg/ml) and DMSO (>50mg/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)