

# 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 G<sub>0</sub>/G<sub>1</sub> transition of the cell cycle (10μM) and causes apoptosis (15μM or greater) in Molt-4 leukemia cells.

Citations: 9

[View Online »](#)

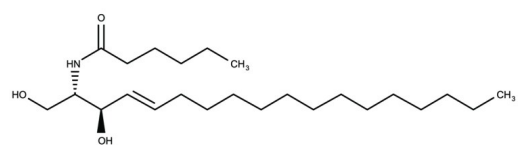
## Ordering Information

[Order Online »](#)

BML-SL110-0005	5mg
BML-SL110-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

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



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-be@enzolifesciences.com](mailto:info-be@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)