

# Sphingosine, D-erythro

## Potent and selective inhibitor of protein kinase C

Common component of cell membranes, often as phosphatidyl conjugates. Is a potent ( $IC_{50}=1-3\text{ }\mu\text{M}$ ) and selective inhibitor of protein kinase C. Inhibition is competitive with diacylglycerol, phorbol dibutyrate and  $Ca^{2+}$  and it also inhibits PKC activation by other lipids. It acts on an equimolar basis with 1,2-dioleoylglycerol and does not affect other kinases such as MLCK and PKA. It is active in intact cells and has been identified as an endogenous constituent in HL-60 cells, neutrophils, rat liver, and brain and mouse tissues. Other activities include inhibition of phosphatidate phosphohydrolase,  $Na^+,K^+$ -ATPase, CTP:phosphocholine cytidyltransferase, calmodulin-dependent enzymes, binding of factor VII to tissue factor, binding of thyrotropin releasing hormone to its receptor and activation of EGF receptor kinase, phospholipase D and casein kinase II.

Citations: 10

[View Online »](#)

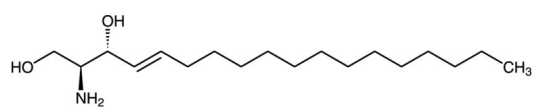
## Ordering Information

[Order Online »](#)

BML-EI155-0025	25mg
BML-EI155-0100	100mg

## 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	D(+)-erythro-1,3-Dihydroxy-2-amino-4- trans-octadecene
Appearance	White to off-white solid.
CAS	123-78-4
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{18}H_{37}NO_2$
MI	14: 8747
MW	299.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (25mg/ml warm) or DMSO (25mg/ml warm).

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)