

Dihydrosphingosine, D-erythro

Sphingosine precursor

Biosynthetic precursor of sphingosine. Inhibitor of protein kinase C (PKC). Blocks phospholipases A₂ (PLA2) and the D-sphingosine precursor.

Citations: 8

[View Online »](#)

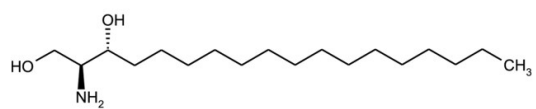
Ordering Information

[Order Online »](#)

BML-SL125-0010	10mg
BML-SL125-0050	50mg

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	D-erythro-Sphingosine, Dihydro-, Sphinganine
Appearance	White to off-white solid.
CAS	764-22-7
Couple Target	Phospholipase, PKC, PLA
Couple Type	Inhibitor
Formula	$C_{18}H_{39}NO_2$
MW	301.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (25mg/ml, warm) or DMSO (25mg/ml, warm).
Technical Info / Product Notes	Note: Product is not sterile.

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)