

# Dihydrosphingosine-1-phosphate, D-erythro

## S1P receptor ligand

Saturated analog of sphingosine-1-phosphate. Induces chemotaxis in HUVECs. May be used as a negative control for intracellular effects of S1P. Rapidly dephosphorylated by a sphingosine phosphatase in cultured human fibroblasts. Ligand for S1P (EDG) receptors.

Citations: 3

[View Online »](#)

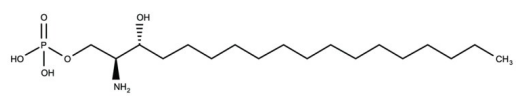
## Ordering Information

[Order Online »](#)

BML-SL143-0001	1mg
----------------	-----

## 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

Appearance	White solid.
CAS	28508-07-8
Couple Target	S1P receptor
Couple Type	Ligand
Formula	$C_{18}H_{40}NO_5P$
MW	381.5
Purity	≥98%
Solubility	Soluble in methanol/THF/water (60/30/10) (0.5mg/ml). Slightly soluble in DMSO, methanol or acetic acid.



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)