

# Sphingomyelin

Ultra-pure

Contains primarily stearic and nervoic acids. Precursor to ceramide second messengers *via* the action of sphingomyelinase.

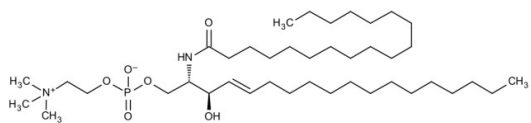
## Ordering Information

[Order Online »](#)

|                |       |
|----------------|-------|
| BML-SL135-0100 | 100mg |
| BML-SL135-0500 | 500mg |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**CAS** 85187-10-6

**Formula**  $C_{41}H_{83}N_2O_6P$

**MW** 731.1 (stearoyl form)

**Purity** ≥98% (TLC)

**Solubility** Soluble in chloroform, ethanol or methanol.

**Source** Purified from bovine spinal cord.

**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)