

3-O-Methyl-Sphingomyelin

Sphingomyelinase inhibitor

3-O-Methyl-Sphingomyelin is a novel selective inhibitor of neutral sphingomyelinase, ($K_i=37\text{ }\mu\text{M}$). It is cell-permeable and does not inhibit acid sphingomyelinase.

Citations: 5

[View Online »](#)

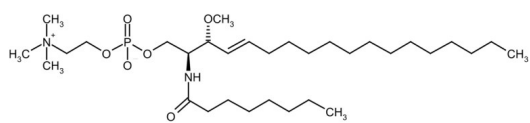
Ordering Information

[Order Online »](#)

BML-SL225-0001	1mg
BML-SL225-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
Couple Target	Sphingomyelinase
Couple Type	Inhibitor
Formula	$C_{32}H_{65}N_2O_6P$
Identity	Determined by NMR.
MW	604.8
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (>25 mg/ml) or DMSO (20 mg/ml@50°C).



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)