

Ganglioside GM1 . sodium salt (bovine brain)

Specific receptor for cholera toxin.

Citations: 3

[View Online »](#)

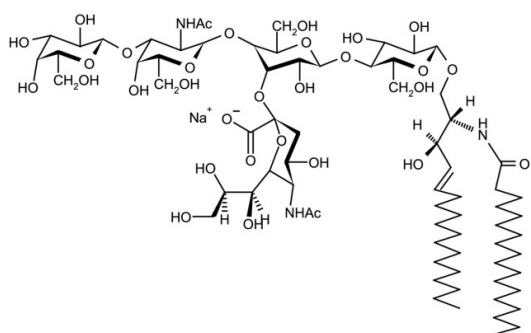
Ordering Information

[Order Online »](#)

ALX-302-001-M001	1mg
ALX-302-001-M005	5mg
ALX-302-001-M010	10mg

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

Alternative Name	GM1 . Na (bovine brain), Monosialoganglioside GM1 . Na (bovine brain)
Appearance	Lyophilized.
CAS	37758-47-7
Couple Target	Cholera toxin
Couple Type	Receptor
Formula	$C_{73}H_{130}N_3O_{31} \cdot Na$
MI	14: 4364
MW	1545.8 . 23.0 (calculated on sphingosine C18:1 and stearic acid)
Purity	≥98% (TLC)
Sequence	[II ³ Neu5AcGgOse ₄ Cer] [β-Gal-(1→3)-β-GalNAc-(1→4)-[α-Neu5Ac-(2→3)]-β-Gal(1→4)-β-Glc-(1→1')-Cer]

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)