

Ganglioside GM3 . sodium salt (bovine brain)

The major ganglioside of the extra-nervous system. Inhibits growth of epidermal cells.

Citations: 3

[View Online »](#)

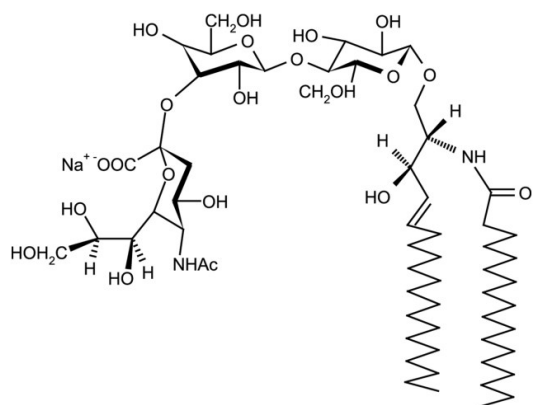
Ordering Information

[Order Online »](#)

ALX-302-005-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GM3 . Na (bovine brain), Monosialoganglioside GM3 . Na (bovine brain)
Appearance	Lyophilized.
CAS	54827-14-4
Formula	$C_{59}H_{107}N_2O_{21} \cdot Na$
MW	1180.5 . 23.0 (calculated on sphingosine C18:1 and stearic acid)
Purity	≥98% (TLC)
Sequence	[II ³ Neu5AcLacCer] [α-Neu5Ac-(2→3)]-β-Gal(1→4)-β-Glc-(1→1')-Cer]
Solubility	Soluble in chlorform/methanol (2:1).

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-
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)