

Ganglioside GD1b . disodium salt (bovine brain)

Citations: 2

[View Online »](#)

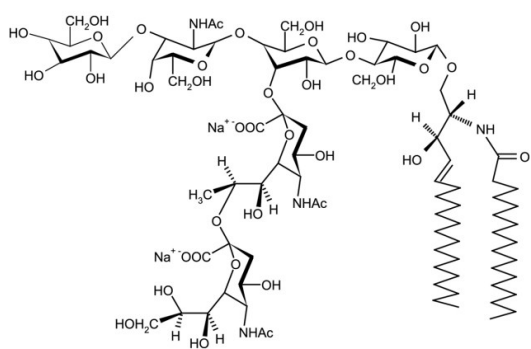
Ordering Information

[Order Online »](#)

ALX-302-009-M001	1mg
------------------	-----

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 Disialoganglioside GD1b . 2Na (bovine brain), GD1b . 2Na (bovine brain)

Appearance Lyophilized.

CAS 19553-76-5

Formula $C_{84}H_{146}N_4O_{39} \cdot 2Na$

MW 1836.1 . 46.0 (calculated on sphingosine C18:1 and stearic acid)

Purity ≥98% (TLC)

Sequence $[II^3(Neu5Ac)_2GgOse_4Cer] [\beta-Gal-(1 \rightarrow 3)-\beta-GalNAc-(1 \rightarrow 4)-[\alpha-Neu5Ac-(2 \rightarrow 8)-\alpha-Neu5Ac-(2 \rightarrow 3)]-\beta-Gal-(1 \rightarrow 4)-\beta-Glc-(1 \rightarrow 1')-Cer]$

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-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)