

Ganglioside GD1a . disodium salt (bovine brain)

PKC inhibitor

Differentiation marker for cell growth. Protein kinase C (PKC) inhibitor. The major ganglioside of the nervous system.

Citations: 5

[View Online »](#)

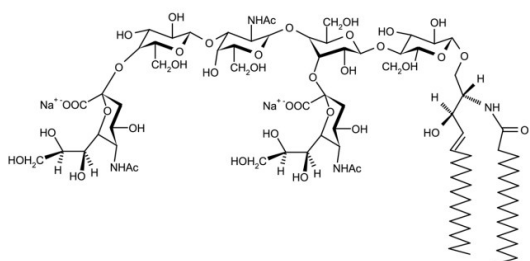
Ordering Information

[Order Online »](#)

ALX-302-007-M001	1mg
ALX-302-007-M005	5mg

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 GD1a . 2Na (bovine brain), GD1a . 2Na (bovine brain)

Appearance Lyophilized.

CAS 12707-58-3

Couple Target PKC

Couple Type Inhibitor

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 $[II^3(Neu5Ac), IV^3(Neu5Ac)GgOse_4Cer] [[\alpha\text{-}Neu5Ac\text{-(}2\rightarrow3\text{)}]\text{-}\beta\text{-Gal}\text{-(}1\rightarrow3\text{)}\text{-}\beta\text{-GalNAc}\text{-(}1\rightarrow4\text{)}\text{-}[\alpha\text{-}Neu5Ac\text{-(}2\rightarrow3\text{)}]\text{-}\beta\text{-Gal}\text{-(}1\rightarrow4\text{)}\text{-}\beta\text{-Glc}\text{-(}1\rightarrow1'\text{)}\text{-Cer}]$

Solubility Soluble in chloroform/methanol/DI water, 2:1:0.1

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@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)