

Asialo-GM1

Citations: 1

[View Online »](#)

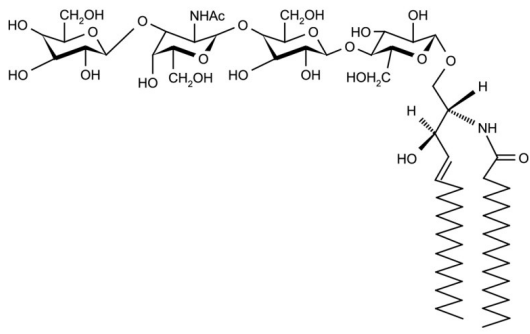
Ordering Information

[Order Online »](#)

ALX-302-013-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	Ceramide tetrahexoside, Gangliotetraosylceramide, Asialoganglioside-GM1
CAS	71012-19-6
Formula	$C_{62}H_{114}N_2O_{23}$
Formulation	Lyophilized.
MW	1255.6 (calculated on sphingosine C18:1 and stearic acid)
Purity	Single Spot (TLC)
Sequence	GgOse ₄ Cer; β-Gal-(1→3)-β-GalNAc-(1→4)-β-Gal-(1→4)-β-Glc-(1→1')-Cer
Source	Semisynthetic (from GM1).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com