

Asialo-GM2

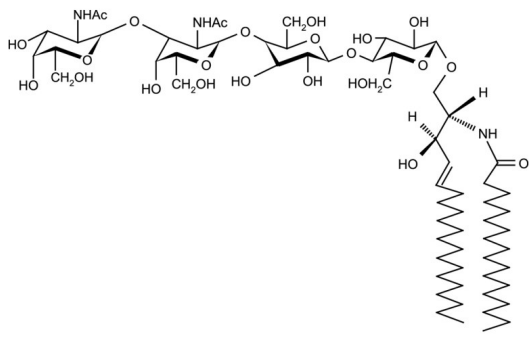
A major neural ganglioside in Tay Sachs and Sandhoff diseases.

Citations: 1

[View Online »](#)

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	Asialoganglioside-GM2, Ceramide trihexoside, Ganglioside GM2
CAS	35960-33-9
Formula	$C_{56}H_{104}N_2O_{18}$
Formulation	Lyophilized.
MW	1093.4 (calculated on sphingosine C18:1 and stearic acid)
Purity	≥98%
Sequence	GgOse ₃ Cer; β-GalNAc-(1→4)-β-Gal-(1→4)-β-Glc-(1→1')-Cer
Source	Semisynthetic (from GM2).

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-fr@enzolifesciences.com

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

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

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