

GM1 Pentasaccharide

. sodium salt

Citations: 5

[View Online »](#)

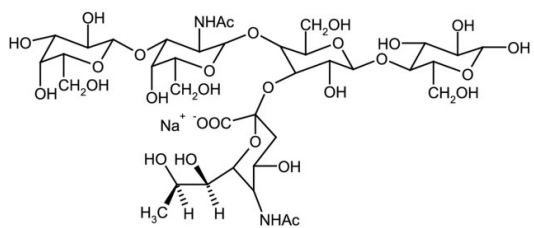
Ordering Information

[Order Online »](#)

ALX-305-006-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

Appearance Lyophilized powder.

CAS 67063-78-9

Formula $C_{37}H_{61}N_2O_{29} \cdot Na$

MW 997.9 . 23.0

Purity ≥95% (1H-NMR)

Sequence $[II^3\text{Neu5AcGgOse}_4] [\beta\text{-Gal-(1}\rightarrow\text{3)}-\beta\text{-GalNAc-(1}\rightarrow\text{4)}-[\alpha\text{-Neu5Ac-(2}\rightarrow\text{3)}]-\beta\text{-Gal-(1}\rightarrow\text{4)}\text{-Glc}]$

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