

Ganglioside GD3 . disodium salt (bovine buttermilk)

Induces mitochondrial permeability transition (MPT) without requiring elevated Ca^{2+} levels and thus triggers Fas-mediated apoptosis.

Citations: 5

[View Online »](#)

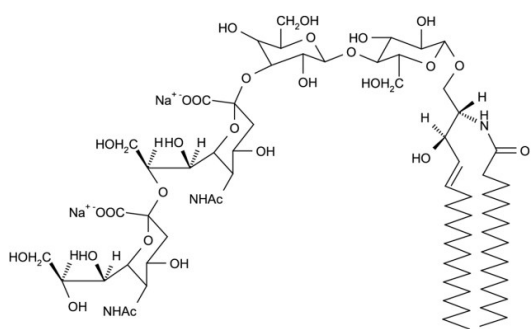
Ordering Information

[Order Online »](#)

ALX-302-010-MC05	0.5mg
ALX-302-010-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 GD3 . 2Na (bovine buttermilk), GD3 . 2Na (bovine buttermilk)

Appearance Lyophilized.

CAS 62010-37-1

Formula $C_{70}H_{123}N_3O_{29} \cdot 2Na$

MW 1470.8 . 46.0 (calculated on sphingosine C18:1 and stearic acid)

Purity ≥98% (TLC)

Sequence $[I\text{I}^3(\text{Neu5Ac})_2\text{LacCer}] [[\alpha\text{-Neu5Ac-(2}\rightarrow\text{8)}\text{-}\alpha\text{-Neu5Ac-(2}\rightarrow\text{3)}]\text{-}\beta\text{-Gal-(1}\rightarrow\text{4)}\text{-}\beta\text{-Glc-(1}\rightarrow\text{1')}\text{-Cer}]$

Solubility Soluble in chloroform/methanol, 2:1

Source Bovine buttermilk

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-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