

# $\alpha$ -Mannosyl Ceramide

## Potent stimulator of NKT cells

Potent stimulator of natural killer T (NKT) cells expressing the invariant V $\alpha$ 19-J $\alpha$ 26 T cell receptor  $\alpha$  chain (V $\alpha$ 19 NKT cells).

Natural killer T cells (NKT cells) are important regulators of the immune system. They share properties of both innate as well as adaptive immunity. They express T cell receptors (TCR) on their surface consisting of an invariable  $\alpha$ -chain and a somewhat variable  $\beta$ -chain. The majority of NKT cells express the invariable  $\alpha$ -chain V $\alpha$ 14-J $\alpha$ 18 (V $\alpha$ 14) in mouse or its homolog in humans V $\alpha$ 24-J $\alpha$ 18 (V $\alpha$ 24). V $\alpha$ 14 NKT cells recognize  $\alpha$ -GalCer presented by CD1d. NKT cells expressing the invariant V $\alpha$ 19-J $\alpha$ 26 TCR  $\alpha$ -chain (V $\alpha$ 19 NKT cells) have been found in peripheral blood lymphocytes and in the liver. They recognize  $\alpha$ -ManCer presented by MR1 (Major histocompatibility complex, class I-related) and respond with production of large amounts of immunoregulatory cytokines.

Citations: 3

[View Online »](#)

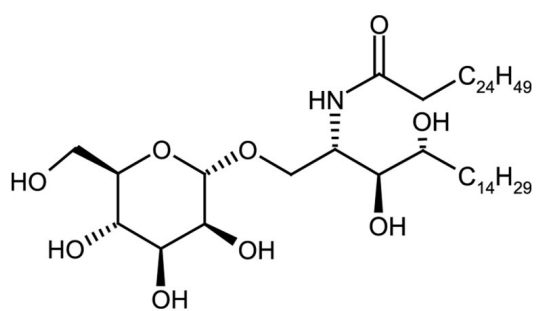
## Ordering Information

[Order Online »](#)

ALX-306-031-M001	1mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After reconstitution, aliquot into glass vials.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	α-ManCer
Appearance	White to brown powder.
Formula	$C_{50}H_{99}NO_9$
MW	858.3
Purity	≥95% (TLC and NMR)
Solubility	Soluble in pyridine or warm DMSO (4mg/ml).



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

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