

N- Butyldeoxynojirimycin . HCl

Inhibitor of glucosylceramide synthetase and glucosidase

A novel inhibitor of glucosylceramide synthase as well as glucosidase I and II. Inhibits glucosylceramide synthase without ceramide accumulation and cell death. It is a potent inhibitor of human immunodeficiency virus (HIV) replication and syncytium formation *in vitro*. Restores chloride current in epithelial CF cells.

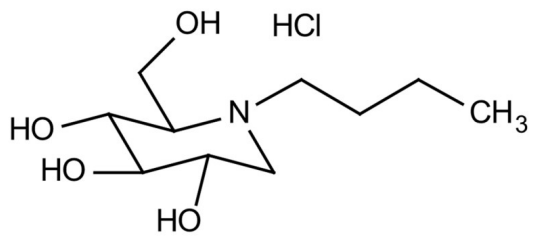
Ordering Information

[Order Online »](#)

BML-SL230-0005	5mg
----------------	-----

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	Miglustat
Appearance	White to off-white solid.
CAS	210110-90-0
Couple Target	Glucosidase, Glucosylceramide synthetase
Couple Type	Inhibitor
Formula	$C_{10}H_{21}NO_4 \cdot HCl$
Identity	Determined by 1H-NMR and MS.
MW	255.7
Purity	≥98% (TLC)
Solubility	Soluble in water.

Last modified: May 29, 2024



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)