

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.

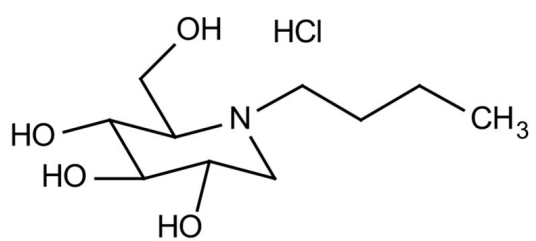
Ordering Information

[Order Online »](#)

BML-SL230-0005	5mg
BML-SL230-0025	25mg

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.



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