

Conduritol B Epoxide

Glycosidase inhibitor

A potent, irreversible inhibitor of various plant β -glucosidases and mammalian glucocerebrosidase. Has been employed in *in vitro* and animal models for Gaucher's disease. It raises intracellular levels of glucosylceramide by inhibiting glucosylceramidase.

Citations: 9

[View Online »](#)

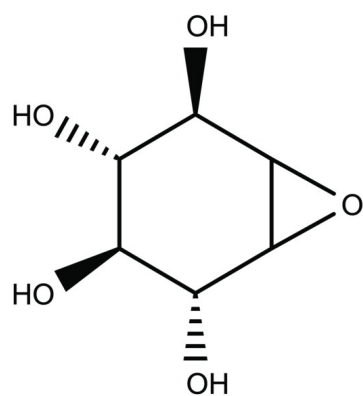
Ordering Information

[Order Online »](#)

BML-S104-0005	5mg
BML-S104-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Product must be stored under argon.
Handling	Hygroscopic. Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	D,L-1,2-Anhydro-myo-inositol
Appearance	White solid.
CAS	6090-95-5
Couple Target	Glucosidase, Glycosidase
Couple Type	Inhibitor
Formula	$C_6H_{10}O_5$
Identity	Determined by ¹ H-NMR and MS.
MW	162.1
Purity	≥96% (TLC)
Solubility	Soluble in water (25mg/ml) or in DMSO.



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