

Swainsonine

Lysosomal alpha-mannosidase inhibitor

Reversibly inhibits the active site of lysosomal α -mannosidase. Also inhibits the growth of tumors and prevents metastasis in murine models. Could inhibit cell proliferation *in vitro* and the growth of human gastric carcinoma *in vivo*. Blocks the processing of high mannose oligosaccharides to form complex type oligosaccharides.[^]Immunomodulator.

Citations: 10

[View Online »](#)

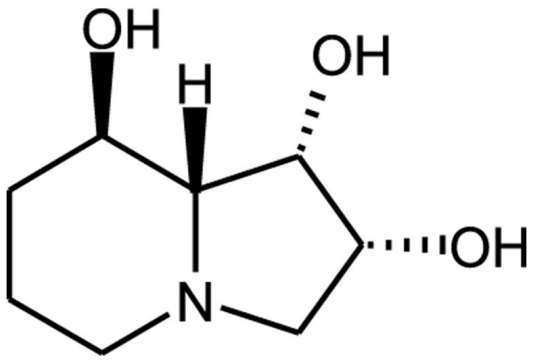
Ordering Information

[Order Online »](#)

BML-S112-0001	1mg
---------------	-----

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 at +4°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	8 α ,8 β -Octahydroindolizidine-1 α ,2 α ,8 β -triol
Appearance	White to off-white solid.
CAS	72741-87-8
Couple Target	Lysosomal alpha-mannosidase, Mannosidase
Couple Type	Inhibitor
Formula	C ₈ H ₁₅ NO ₃
MW	173.2
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, DMSO, methanol or water (10mg/ml).

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-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