

# Swainsonine

## Lysosomal alpha-mannosidase inhibitor

Reversibly inhibits the active site of lysosomal  $\alpha$ -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.<sup>^</sup>Immunomodulator.

Citations: 10

[View Online »](#)

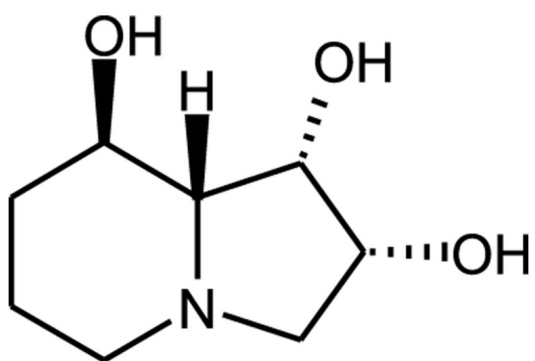
## Ordering Information

[Order Online »](#)

BML-S112-0001	1mg
BML-S112-0005	5mg

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