

# Castanospermine

## Glucosidase inhibitor

Plant alkaloid. Potent inhibitor of  $\alpha$ - and  $\beta$ -glucosidases. Exhibits antiviral properties. Inhibits thyroglobulin secretion. Antitumor activity. Inhibits protein glycosylation. Prevents angiogenesis.

Citations: 15

[View Online »](#)

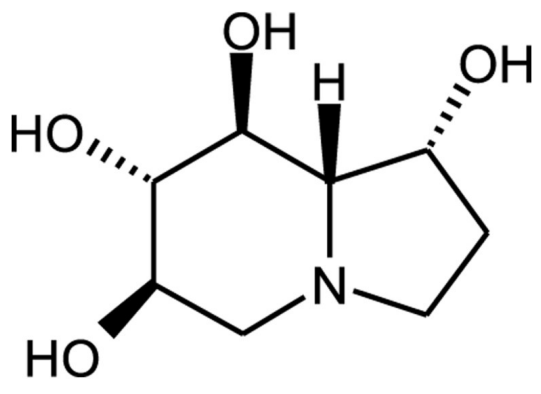
## Ordering Information

[Order Online »](#)

BML-S107-0100	100mg
BML-S107-0500	500mg

## 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

Appearance	White to off-white crystalline solid.
CAS	79831-76-8
Couple Target	Glucosidase, Glycosidase
Couple Type	Inhibitor
Formula	$C_8H_{15}NO_4$
MI	14: 1896
MW	189.2
Purity	≥95% (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-  
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)