

# Manzamine A

GSK-3 inhibitor

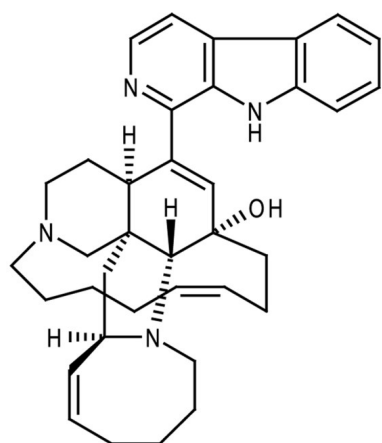
$\beta$ -Carboline alkaloid with anti-inflammatory, antimalarial, antifungal, anti-HIV-1 and insecticidal activity. Also active against the Gram-positive bacteria *Bacillus subtilis* and *Staphylococcus aureus*. Inhibits GSK-3. Has native fluorescence, highest when solubilized in methanol (Ex: 340nm, Em: 387nm).

Citations: 8

[View Online »](#)

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

**Appearance** Off white to yellow solid.

**CAS** 104196-68-1

**Couple Target** GSK

**Couple Type** Inhibitor

**Formula**  $C_{36}H_{44}N_4O$

**MW** 548.8

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO, methanol or 100% ethanol.

**Source** Isolated from *Xestospongia* sp.

**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-  
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)