

# Veratridine

Na<sup>+</sup> channel activator

A Na<sup>+</sup> channel activator. Does not affect Na<sup>+</sup>-Ca<sup>2+</sup> exchange. Veratridine acts at neurotoxin receptor site 2 and preferentially binds to activated Na<sup>+</sup> channels causing persistent activation. It prevents inactivation and shifts activation to more negative potentials. Induces cell death in bovine chromaffin cells.

Citations: 16

[View Online »](#)

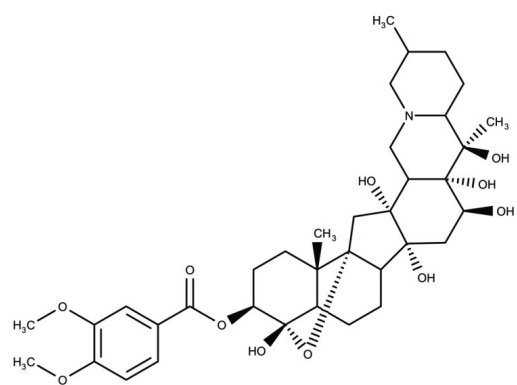
## Ordering Information

[Order Online »](#)

BML-NA125-0010	10mg
----------------	------

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

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White to yellow solid.
CAS	71-62-5
Couple Target	Sodium channel
Couple Type	Activator
Formula	$C_{36}H_{51}NO_{11}$
MW	673.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)