

# $\alpha$ -Latrotoxin

## Neurotoxin

Causes massive neurotransmitter release from a wide variety of central and peripheral synaptic junctions of vertebrates using  $\text{Ca}^{2+}$ -dependent and  $\text{Ca}^{2+}$ -independent pathways. A useful pharmacological tool in the studies of synaptic vesicles exocytosis of different neurotransmitters.

Citations: 13

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-630-027-C040	40µg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution is stable for 1 week when stored at +4°C or for at least one year at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

CAS	65988-34-3
Formulation	Lyophilized.
MW	~130kDa.
Purity	≥97%
Purity Detail	Migrates as one distinct band when run on a 5-25% gradient polyacrylamide gel according to the method of Laemmli (Nature 227, 680 (1970)).
Quality Control	Bioassay: Test of the ability to stimulate neurotransmitter release in both Ca <sup>2+</sup> -free and Ca <sup>2+</sup> -containing media according to Valtorta. Effective concentration: 100pM-1nM.
Quantity	Determined by Lowry and Pierce method.
RTECS	OE9020000
Reconstitution	Reconstitute in 0.5ml of distilled cold water. Stir very gently. Do not shake or vortex. Add 0.5 ml glycerol. Stock solution received contains 50% glycerol and 300nM α-latrotoxin.
Source	Isolated from <i>Latrodectus tredecimguttatus</i> .
Technical Info / Product Notes	<b>Attention:</b> For best results, only proceed according to the given instructions. Dissolution and/or storage of this toxin contrary to these instructions may affect the stability and biological activity of this compound.



ENZO LIFE SCIENCES  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**P23631**  
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)