

# Convulxin

## Heterodimeric C-type lectin

Heterodimeric C-type lectin. Activates mammalian platelets via binding and clustering of p62/GPVI-receptors under physiological conditions.

Citations: 28

[View Online »](#)

### Ordering Information

[Order Online »](#)

ALX-350-100-C050	50µg
------------------	------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for 8 hours at +20°C, for 2 days at +4°C or for 1 month when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Activity	<50ng/ml (minimal concentration to induce maximum activation of washed human platelets).
Application Notes	Can be used for studies on platelet receptors.
CAS	37206-04-5
Formulation	Lyophilized.
MW	~84kDa.
Purity	≥90% (SDS-PAGE)
Purity Detail	Partially purified by multi-step chromatography.
Reconstitution	Reconstitute in 10mM HEPES buffer, pH 7.2. (We recommend adding a carrier protein such as BSA to the reconstitution buffer and any subsequent dilution buffers to prevent absorption on plastic surfaces.)
Solubility	Soluble in organic buffer (HEPES).
Source	Isolated from <i>Crotalus durissus terrificus</i> snake venom.

## UniProt ID

O93426



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)

Last modified: May 29, 2024