

# Chlorotoxin

Cl<sup>-</sup> channel blocker

Potent neurotoxin. Small conductance Cl<sup>-</sup> channel blocker. Highly specific marker for glioma cells and tumors of neuroectodermal origin. Inhibits the activity of matrix metalloproteinase-2 (MMP-2).

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-630-069-C500	500µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Aliquots are stable for a few weeks when stored at -20°C.
<b>Handling</b>	After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	Lyophilized.
<b>CAS</b>	163515-35-3
<b>Formula</b>	$C_{158}H_{249}N_{53}O_{47}S_{11}$
<b>MW</b>	3995.8
<b>Purity</b>	≥98% (TLC, HPLC)
<b>Quantity</b>	Vial contains exactly 0.55mg (0.14μmol)
<b>Reconstitution</b>	Reconstitute by injecting 1.4ml of distilled water into the vial using a calibrated syringe. Then dissolve the contents thoroughly to receive a 10-4 M solution.
<b>Sequence</b>	H-Met-Cys-Met-Pro-Cys-Phe-Thr-Thr-Asp-His-Gln-Met-Ala-Arg-Lys-Cys-Asp-Asp-Cys-Cys-Gly-Gly-Lys-Gly-Arg-Gly-Lys-Cys-Tyr-Gly-Pro-Gln-Cys-Leu-Cys-Arg-NH <sub>2</sub> (Disulfide bonds between Cys <sup>2</sup> -Cys <sup>19</sup> , Cys <sup>5</sup> -Cys <sup>28</sup> , Cys <sup>16</sup> -Cys <sup>33</sup> and Cys <sup>20</sup> -Cys <sup>35</sup> )
<b>Solubility</b>	Soluble in water
<b>Source</b>	Synthetic.



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)