

# Dendrotoxin-I (Dendroaspis p. polylepis), (purified)

## Ultra-pure

K<sup>+</sup> channel blocker. Blocks K<sub>V</sub>1.2 and K<sub>V</sub>1.1 channels (IC<sub>50</sub>=0.13 and 3.1nM in oocytes respectively, with higher values for mammalian cells), as well as heteromultimeric channels comprised of K<sub>V</sub>1.2 and K<sub>V</sub>1.1 with other K<sub>V</sub>1 isoforms.

**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	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

CAS	107950-33-4
Formulation	Lyophilized.
MW	7149
Purity	≥98% (HPLC)
Solubility	Soluble in aqueous buffers (pH 7.5).
UniProt ID	P00979

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)