

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

Ultra-pure

K⁺ channel blocker. Blocks K_v1.2 and K_v1.1 channels (IC₅₀=0.13 and 3.1nM in oocytes respectively, with higher values for mammalian cells), as well as heteromultimeric channels comprised of K_v1.2 and K_v1.1 with other K_v1 isoforms.

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| BML-KC163-0140 | 140µg |
|----------------|-------|

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com