

α -Dendrotoxin

K⁺ channel blocker

α -Dendrotoxin is isolated from *Dendroaspis angusticeps* snake venom, by modification of the procedures of Harvey and Benishin and purified to homogeneity. α -Dendrotoxin blocks K_V1.1 and K_V1.2 channels (IC₅₀ = 0.4 to 4 and 1.1 to 12 nM in oocytes respectively, with higher values for mammalian cells). In addition, the toxin was shown to block also K_V1.6 (IC₅₀ = 9-25 nM).

Ordering Information

[Order Online »](#)

BML-KC167-0140	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

Appearance Lyophilized solid.

CAS 74504-53-3

MW 7048.1

Purity ≥98% (HPLC)

Sequence Gln-Pro-Arg-Arg-Lys-Leu-Cys-Ile-Leu-His-Arg-Asn-Pro-Gly-Arg-Cys-Tyr-Asp-Lys-Ile-Pro-Ala-Phe-Tyr-Tyr-Asn-Gln-Lys-Lys-Lys-Gln-Cys-Glu-Arg-Phe-Asp-Trp-Ser-Gly-Cys-Gly-Gly-Asn-Ser-Asn-Arg-Phe-Lys-Thr-Ile-Glu-Glu-Cys-Arg-Arg-Thr-Cys-Ile-Gly

Solubility Soluble in any aqueous buffer. Please centrifuge all product preparations at 10000 x g for 5 minutes before reconstitution.



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)