

Apamin

K⁺ channel blocker

Originally isolated from *Apis mellifera*. A small conductance Ca²⁺-activated K⁺ channel blocker.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-151-013-MC05	0.5mg
------------------	-------

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Lyophilized.

CAS 24345-16-2

Formula $C_{79}H_{131}N_{31}O_{24}S_4$

MI **14:** 725

MW 2027.3

Purity ≥95% (HPLC)

RTECS CD6899900

Sequence H-Cys-Asn-Cys-Lys-Ala-Pro-Glu-Thr-Ala-Leu-Cys-Ala-Arg-Arg-Cys-Gln-Gln-His-NH₂
(Disulfide bonds between Cys¹-Cys¹¹ and Cys³-Cys¹⁵)

Solubility Soluble in 0.05M acetic acid (5mg/ml); slightly soluble in water (1mg/ml).

Source Synthetic.

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)