

Verruculogen

Cell cycle inhibitor

Neurotoxin that acts as an inhibitor of the M phase of the mammalian cell cycle. Inhibits Ca^{2+} -activated K^+ channels. Decreases GABA levels in CNS.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-380-056-M001	1mg
ALX-380-056-M005	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 White to off-white powder.

CAS 12771-72-1

Couple Target Potassium channel

Couple Type Inhibitor

Formula $C_{27}H_{33}N_3O_7$

MW 511.6

Purity ≥95% (HPLC, TLC)

RTECS JH4920000

Solubility Soluble in ethyl acetate, DMSO, 100% ethanol or methanol.

Source Isolated from *Penicillium verruculosum*.

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)