

Penitrem A

K⁺ channel blocker

Fungal mycotoxin. Selective, irreversible blocker of the smooth muscle high conductance Ca²⁺-activated K⁺ (maxi-K) channel (100% block at 10 nM). Increases spontaneous release of GABA and aspartate from cerebrocortical synaptosomes in rat neuromuscular junction.

Citations: 7

[View Online »](#)

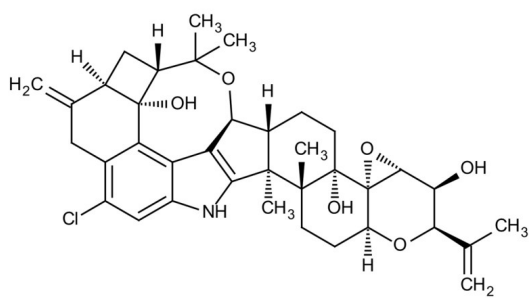
Ordering Information

[Order Online »](#)

BML-KC157-0001	1mg
BML-KC157-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NSC 354845, Tremortin A
Appearance	White solid.
CAS	12627-35-9
Couple Target	Potassium channel
Couple Type	Blocker
Formula	$C_{37}H_{44}ClNO_6$
MW	634.2
Purity	≥95% (HPLC, TLC)
RTECS	RY7535000
Solubility	Soluble in organic solvents such as methanol, acetone or DMSO.

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-
fr@enzolifesciences.com](mailto:info-fr@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)