

# Penitrem A

K<sup>+</sup> channel blocker

Fungal mycotoxin. Selective, irreversible blocker of the smooth muscle high conductance Ca<sup>2+</sup>-activated K<sup>+</sup> (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 »](#)

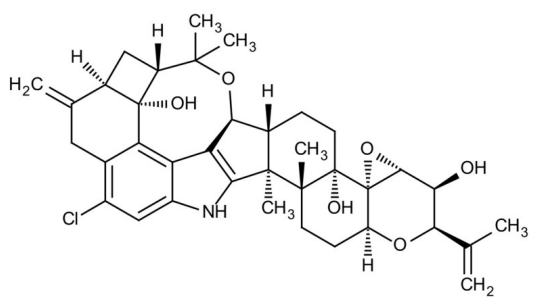
## Ordering Information

[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.



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)