

Paxilline

K⁺ channel blocker

Paxilline is an indole alkaloid isolated from *Penicillium paxilli*. It is a selective and reversible blocker of the smooth muscle high conductance Ca²⁺-activated K⁺ (maxi-K) channel (70% block at 10 nM). In neutrophils, Paxilline blockage of BK_{Ca} channels decreased the alkalinization of phagocytic vacuoles and inhibited the microbicidal process. Inhibitor of the cerebellar inositol 1,4,5-triphosphate (InsP(3)) receptor.

Citations: 9

[View Online »](#)

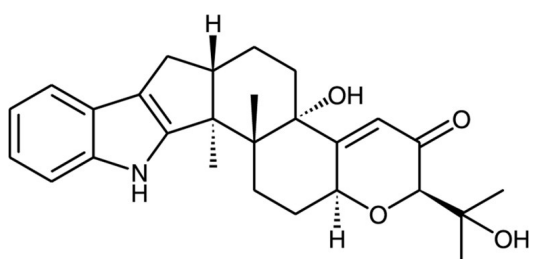
Ordering Information

[Order Online »](#)

BML-KC155-0005	5mg
BML-KC155-0025	25mg

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

Appearance	White solid.
CAS	57186-25-1
Couple Target	IP3 receptor, Potassium channel
Couple Type	Blocker, Ligand
Formula	$C_{27}H_{33}NO_4$
MW	435.6
Purity	≥97% (HPLC)
RTECS	DJ2830000
Solubility	Soluble in DMSO (50mg/ml).



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)