

Isotetrandrine

Phospholipase inhibitor

Isotetrandrine is a biscoclaurine alkaloid which inhibits G-protein mediated activation of phospholipase A₂ (PLA₂) but not phospholipase C (PLC) or D (PLD). Cell permeable.

Citations: 5

[View Online »](#)

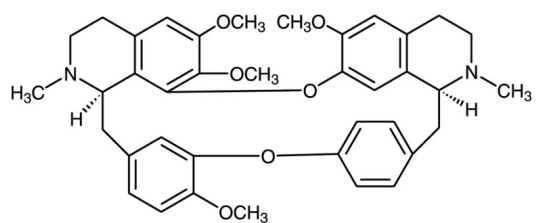
Ordering Information

[Order Online »](#)

BML-G520-0001	1mg
---------------	-----

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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Berbamine methyl ether, 6,6',7,12-Tetramethoxy-2,2'-dimethyl-berbaman
Appearance	Off-white solid.
CAS	477-57-6
Couple Target	Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{38}H_{42}N_2O_6$
MW	622.7
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (20mg/ml) or ethanol (25mg/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)