Isotetrandrine

Phospholipase inhibitor

Isotetrandrine is a biscoclaurine alkaloid which inhibits G-protein mediated activation of phospholipase A_2 (PLA₂) but not phospholipase C (PLC) or D (PLD). Cell permeable.

Citations: 5

View Online »

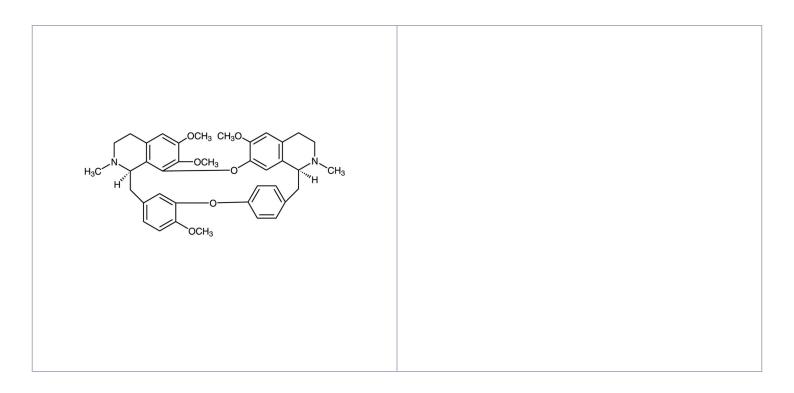
Ordering Information

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)

Soluble in DMSO (20mg/ml) or ethanol (25mg/ml).

Last modified: May 29, 2024

