

Vinpocetine

PDE inhibitor

Selective inhibitor of Ca^{2+} /calmodulin-dependent phosphodiesterase (PDE I, $\text{IC}_{50}=20\text{ }\mu\text{M}$).

Citations: 5

[View Online »](#)

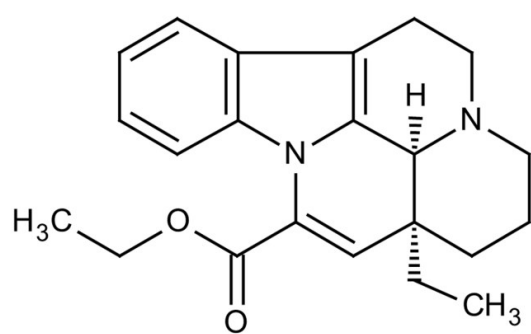
Ordering Information

[Order Online »](#)

BML-PD185-0020	20mg
BML-PD185-0100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White crystalline solid.

CAS 42971-09-5

Couple Target PDE

Couple Type Inhibitor

Formula $C_{22}H_{26}N_2O_2$

Identity Determined by IR.

MI 14: 9991

MW 350.5

Purity ≥98% (Assay)

RTECS JW4792000

Solubility Soluble in DMSO or acetone.

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)