

Trequinsin

PDE inhibitor

Ultrapotent, highly selective inhibitor of cGMP-inhibited phosphodiesterase (PDE III, $IC_{50}=0.3nM$) *in vitro*. Potentiates adenosine-stimulated cAMP accumulation.

Citations: 5

[View Online »](#)

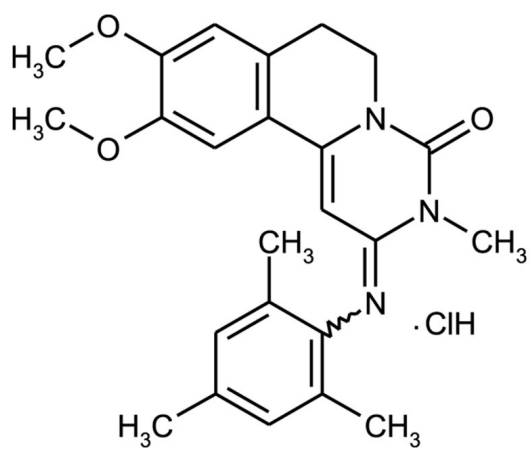
Ordering Information

[Order Online »](#)

BML-PD180-0010	10mg
----------------	------

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 at +20°C. Solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from moisture.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	9,10-Dimethoxy-2-mesitylimino-3-methyl-2,3,6,7-tetrahydro-4H-pyrimido-[6,1a]-isoquinolin-4-one . HCl, HL 725
Appearance	White to pale yellow solid.
CAS	78416-81-6
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_{24}H_{27}N_3O_3 \cdot HCl$
Identity	Determined by 1H-NMR and LC-MS.
MW	405.5 . 36.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (20mg/ml); slightly soluble in water.

Technical Info / Product Notes

Note: Product is not sterile.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024