

8-Methoxymethyl- IBMX

PDE inhibitor

Selective inhibitor of Ca^{2+} -calmodulin-dependent phosphodiesterase (PDE I, $\text{IC}_{50}=4\text{ }\mu\text{M}$).

Citations: 2

[View Online »](#)

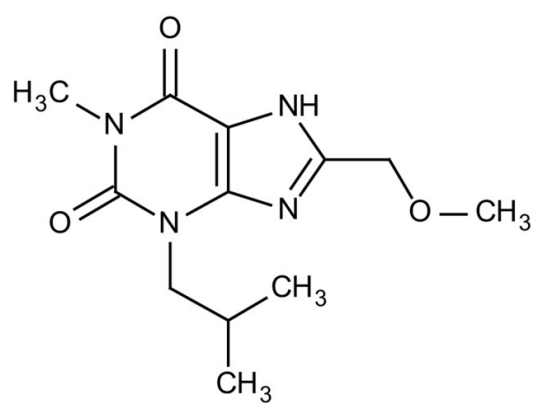
Ordering Information

[Order Online »](#)

BML-PD150-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Use, as supplied, at room temperature for up to 2 years. Store dilutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	78033-08-6
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_{12}H_{18}N_4O_3$
Identity	Determined by 1H -NMR and MS.
MW	266.3
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (5mg/ml) or DMSO (5mg/ml)

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)