

Cilostamide

PDE inhibitor

Phosphodiesterase PDE3 inhibitor ($IC_{50}=5$ nM). Blocks PKB/Akt signaling pathway by inhibition of Akt-activated PDE3B. Reverses the established effects of leptin on food intake and body weight. Blocks the leptin-induced tyrosine phosphorylation of STAT3 at the hypothalamic level.

Citations: 9

[View Online »](#)

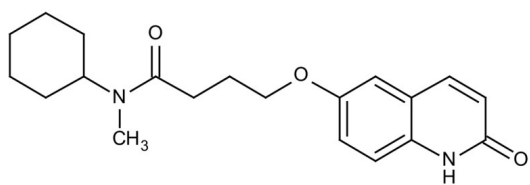
Ordering Information

[Order Online »](#)

BML-PD125-0005	5mg
BML-PD125-0025	25mg

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 room temperature. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light and oxygen. Packaged under inert gas.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	OCP 3689, OPS 3689, N-Cyclohexyl-N-methyl-4-(1,2-dihydro-2-oxo-6-quinolyloxy)butyramide
Appearance	White to off-white solid.
CAS	68550-75-4
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_{20}H_{26}N_2O_3$
Identity	Determined by NMR.
MW	342.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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