

Milrinone

PDE3 inhibitor

Selective inhibitor of phosphodiesterase 3 (PDE3, $IC_{50}=0.3\mu M$).

Citations: 8

[View Online »](#)

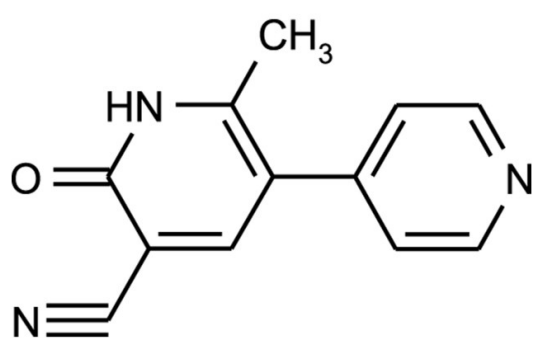
Ordering Information

[Order Online »](#)

ALX-270-083-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,6-Dihydro-2-methyl-6-oxo-3,4'-bipyridine-5-carbonitrile

Appearance White to light red solid.

CAS 78415-72-2

Couple Target PDE

Couple Type Inhibitor

Formula $C_{12}H_9N_3O$

MI 14: 6197

MW 211.2

Purity ≥97% (HPLC)

RTECS DW1762000

Solubility Soluble in DMSO (20mg/ml) or warm 100% ethanol (3mg/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)