

PJ-34

PARP inhibitor

Potent, water soluble poly(ADP-ribose) polymerase (PARP) inhibitor (EC_{50} =20nM compared to EC_{50} =200 μ M of the prototypical PARP inhibitor 3-aminobenzamide (Prod. No. ALX-270-044)). Inhibits peroxynitrite (Prod. No. ALX-400-036)-induced cell necrosis (EC_{50} =20nM). Has significant, dose-dependent, anti-inflammatory effects in a variety of local inflammation models and provides cardioprotection by decreasing myocardial infarct size.

Citations: 32

[View Online »](#)

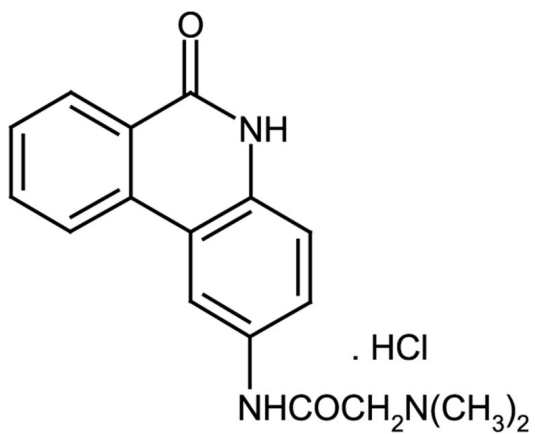
Ordering Information

[Order Online »](#)

ALX-270-289-M001	1mg
ALX-270-289-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2 years after receipt when stored at +4C° (desiccated). Stock solutions are stable for 3 weeks when stored at room temperature.
Handling	Hygroscopic.
Short Term Storage	Ambient
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(6-Oxo-5,6-dihydro-phenanthridin-2-yl)-N,N-dimethylacetamide . HCl
Appearance	White to light brown or yellow powder.
CAS	344458-15-7
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_{17}H_{17}N_3O_2 \cdot HCl$
Identity	Determined by 1H-NMR.
MW	295.2 . 36.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 30mg/ml) or water (up to 20mg/ml).
Source	Synthetic.

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-
be@enzolifesciences.com](mailto:info-be@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)