

# EB-47 . dihydrochloride . dihydrate

## PARP-1 inhibitor

Very potent and water soluble PARP-1 inhibitor ( $IC_{50}$ =45nM, 100% inhibition at 200nM). Shows cytoprotective effects against oxidative damage in cells and in *in vivo* models of reperfusion injury and inflammation.

Citations: 3

[View Online »](#)

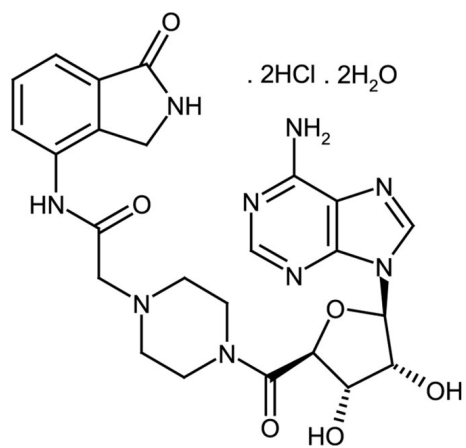
## Ordering Information

[Order Online »](#)

ALX-270-383-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	1-Piperazineacetamide,4-[1-(6-amino-9H-purin-9-yl)-1-deoxy-β-D-ribofuranuron]-N-(2,3-dihydro-1H-isoindol-4-yl)-1-one . 2HCl . 2H <sub>2</sub> O
Appearance	White solid.
Couple Target	PARP
Couple Type	Inhibitor
Formula	C <sub>24</sub> H <sub>27</sub> N <sub>9</sub> O <sub>6</sub> . 2HCl . 2H <sub>2</sub> O
Identity	Determined by 1H-NMR.
MW	537.5 . 73.0 . 36.0
Purity	≥95% (HPLC)
Solubility	Soluble in water or DMSO.

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)