

# ABT-888

## PARP inhibitor

Potent inhibitor of PARP-1 and PARP-2 (potency  $\leq 5\text{nM}$  *in vitro*). Does not inhibit other NAD-binding enzymes. Has minimal CYP450 inhibition and induction. Shows broad spectrum of chemo- and radiopotentialiation. Is toxic to both oxic and hypoxic cells. Enantiomeric purity  $\geq 97\%$  suitable for *in vivo* studies. Does not show inherent cytotoxicity and shows no single agent activity in tumor models. Has excellent bioavailability and good blood-brain permeation. Increases tumor growth delay resulting from radiation and DNA-damaging agents.

Citations: 34

[View Online »](#)

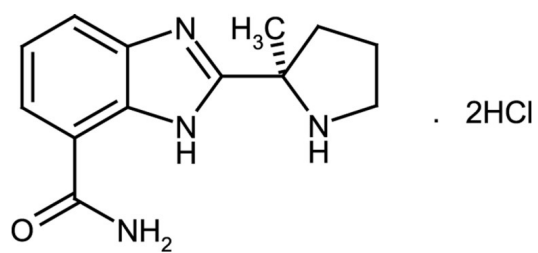
## Ordering Information

[Order Online »](#)

ALX-270-444-M001	1mg
ALX-270-444-M005	5mg
ALX-270-444-M025	25mg

## 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** 2-[(2R)-2-Methylpyrrolidin-2-yl]-1H-benzimidazole-4-carboxamide . dihydrochloride

**Appearance** Colorless to white crystalline solid.

**CAS** 912445-05-7

**Couple Target** PARP

**Couple Type** Inhibitor

**Formula**  $C_{13}H_{16}N_4O \cdot 2HCl$

**Identity** Identity determined by <sup>1</sup>H-NMR.

**MW** 244.3 . 73.0

**Purity** ≥98% (HPLC)

**Solubility** Soluble in water or DMSO.



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)