

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 radiopotentiality. 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 »](#)

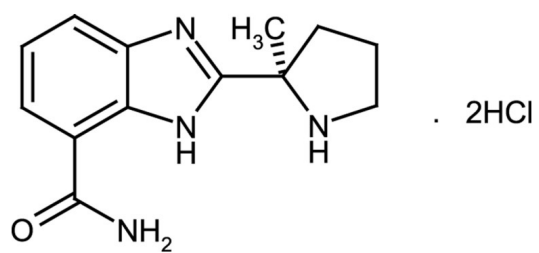
Ordering Information

[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 1H-NMR.

MW 244.3 . 73.0

Purity ≥98% (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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)