

4-Amino-1,8-naphthalimide

PARP inhibitor

Potent inhibitor of poly (ADP-ribose) polymerase-1 (PARP-1) (IC_{50} = 0.18 μ M). Reduces ischemia-reperfusion injury in the heart and skeletal muscle.

Citations: 9

[View Online »](#)

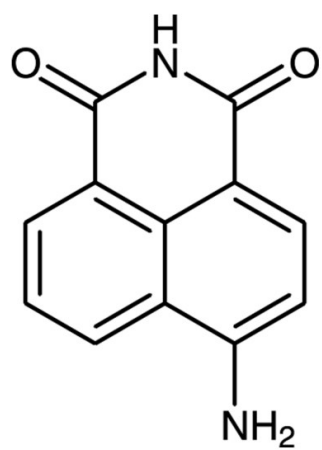
Ordering Information

[Order Online »](#)

BML-AP101-0020	20mg
BML-AP101-0100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-ANI
Appearance	Orange solid.
CAS	1742-95-6
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_{12}H_8N_2O_2$
MW	212.2
Purity	≥96% (NMR)
Solubility	Soluble in DMSO (5mg/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)