

D PQ

PARP-1 inhibitor

Very potent poly(ADP-ribose) polymerase-1 (PARP-1) inhibitor.

Citations: 13

[View Online »](#)

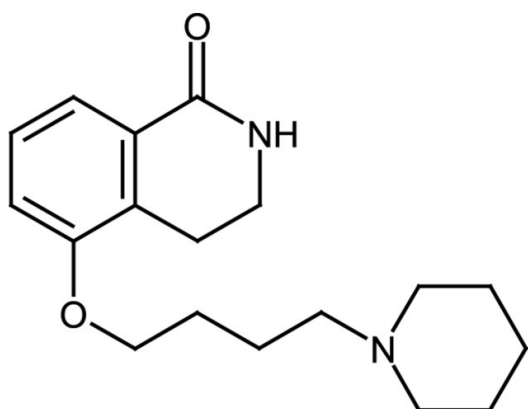
Ordering Information

[Order Online »](#)

ALX-270-221-M001	1mg
ALX-270-221-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for up to 1 month when stored at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white to light brown solid.
CAS	129075-73-6
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_{18}H_{26}N_2O_2$
Identity	Determined by ¹ H-NMR.
MW	302.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO; slightly soluble in 100% ethanol; insoluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com