

UPF1035

PARP-2 inhibitor

PARP-2 specific inhibitor with 60-fold selectivity for PARP-2 ($IC_{50}=0.15 \pm 0.04\mu M$) over PARP-1 ($IC_{50}=9.0 \pm 0.7\mu M$). Can be used for further characterization of PARP-2 in pathophysiological conditions.

Citations: 2

[View Online »](#)

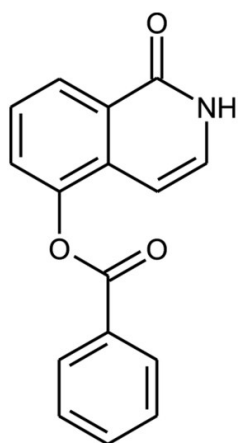
Ordering Information

[Order Online »](#)

ALX-270-498-M005	5mg
------------------	-----

Manuals, SDS & CoFA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Not stable in basic conditions.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Benzoyloxy-3,4-dihydroisoquinolin-1(2H)-one
CAS	370872-09-6
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_{16}H_{11}NO_3$
MW	265.3
Purity	≥95% (HPLC)
Solubility	Soluble in 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-
be@enzolifesciences.com](mailto:info-be@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)