

TIQ-A

PARP-1 inhibitor

Potent poly(ADP-ribose) polymerase-1 (PARP-1) inhibitor (IC_{50} =450nM).
Neuroprotectant.

Citations: 2

[View Online »](#)

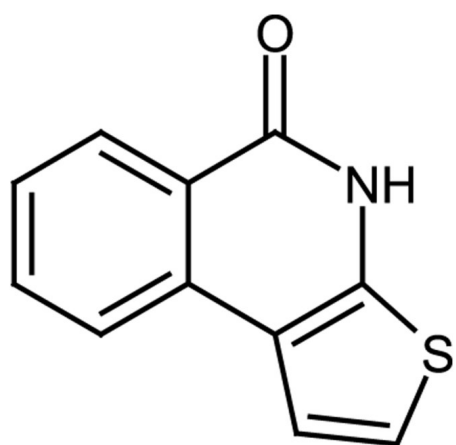
Ordering Information

[Order Online »](#)

ALX-270-365-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Thieno[2,3-c]isoquinolin-5-one

CAS 420849-22-5

Couple Target PARP

Couple Type Inhibitor

Formula C₁₁H₇NOS

MW 201.2

Purity ≥96% (HPLC)

Solubility Soluble in DMSO.

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-be@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