

DR2313

PARP inhibitor

Potent, water soluble competitive PARP inhibitor (IC_{50} =0.20 μ M and 0.24 μ M for PARP-1 and PARP-2 respectively).

Citations: 1

[View Online »](#)

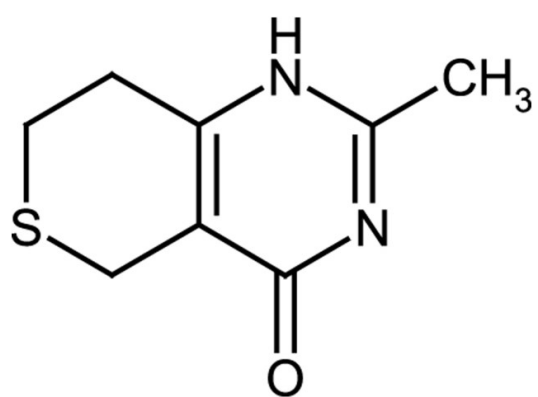
Ordering Information

[Order Online »](#)

ALX-270-452-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Methyl-3,5,7,8-tetrahydrothiopyrano[4,3-d]pyrimidine-4-one
Appearance	White to off-white crystalline solid.
CAS	284028-90-6
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_8H_{10}N_2OS$
MW	182.2
Purity	≥98% (HPLC (UV))
Solubility	Soluble in DMSO, methanol or water.

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)