

NU1025

PARP-1 inhibitor

A potent poly(ADP-ribose) polymerase 1 (PARP-1) inhibitor ($IC_{50}=400nM$) that potentiates the cytotoxicity of various DNA-active agents, including the DNA-methylating compound MTIC, the DNA strand break-inducing drug temozolomide (Prod. No. ALX-420-044), topotecan (Prod. No. ALX-350-133), bleomycin (Prod. No. BML-AP302), and ionizing radiation effective in murine L1210 leukemia cells, CHO, and in a variety of human tumor cell lines.

Citations: 11

[View Online »](#)

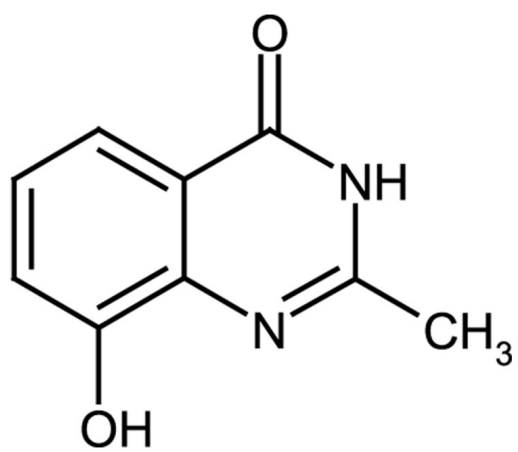
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-270-370-M005 | 5mg |
| ALX-270-370-M025 | 25mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. |
| Handling | Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C. |
| Short Term Storage | -20°C |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|--|
| Alternative Name | 8-Hydroxy-2-methylquinazoline-4-one |
| Appearance | Off-white to purple solid. |
| CAS | 90417-38-2 |
| Couple Target | PARP |
| Couple Type | Inhibitor |
| Formula | $C_9H_8N_2O_2$ |
| MW | 176.2 |
| Purity | ≥98% (HPLC) |
| Reconstitution | Reconstitute with DMSO. |
| Solubility | Soluble in DMSO (25mg/ml). This solution can be further diluted 1:100 in tissue culture medium before use. |

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