

ICRF-193

Topoisomerase inhibitor

A topoisomerase II inhibitor that does not cause DNA damage. Induces apoptosis in several cell lines including K562 and Molt-4 cells. Induces a G2 checkpoint that is associated with an ATR-dependent inhibition of polo-like kinase 1 (plk1) activity and a decrease in cyclin B1 phosphorylation.

Citations: 8

[View Online »](#)

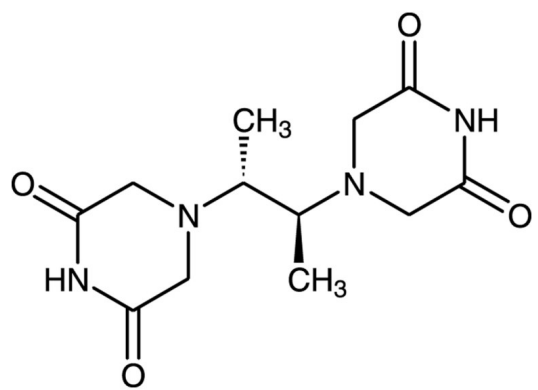
Ordering Information

[Order Online »](#)

BML-GR332-0001	1mg
BML-GR332-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Meso-4,4'-(3,2-butanediyl)-bis(2,6-piperazinedione)

Appearance Off-white to pale yellow solid.

CAS 21416-88-6

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{12}H_{18}N_4O_4$

MW 282.3

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (4mg/ml).

Last modified: June 16, 2025



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)