

# 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 »](#)

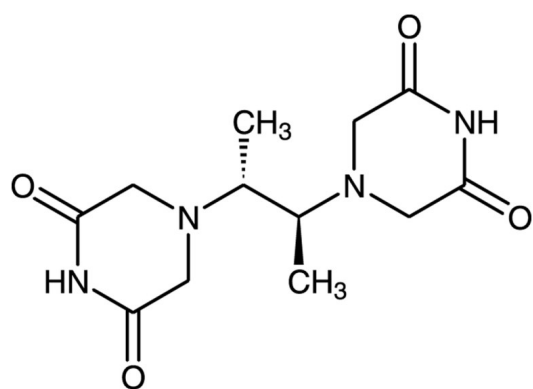
## Ordering Information

[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).



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)