

# PRIMA-1

## p53 activator

Novel cell-permeable p53 reactivating agent. Restores sequence-specific DNA binding and the active conformation to mutant p53 (in 13 of 14 p53 mutants tested) *in vitro* and in living cells. Induces apoptosis in human tumor cell lines and displays antitumor activity in mice.

Citations: 1

[View Online »](#)

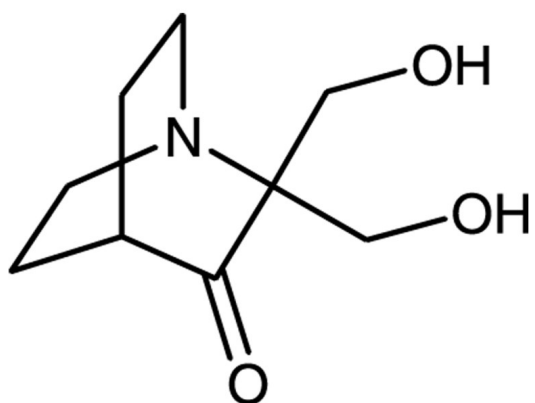
## Ordering Information

[Order Online »](#)

BML-GR333-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store stock solutions at -20°C for up to 3 months. Storage of aqueous solutions is not recommended.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2,2-bis(Hydroxymethyl)-1-azabicyclo[2.2.2]octan-3-one, 2,2-bis(Hydroxymethyl)-3-quinuclidinone
Appearance	White to off-white solid.
CAS	5608-24-2
Couple Target	p53
Couple Type	Activator
Formula	$C_9H_{15}NO_3$
MW	185.2
Purity	≥98% (TLC)
Solubility	Soluble at 25 mg/ml in ethanol, 100% DMSO, and DMF. Soluble at 10 mg/ml in water.
Technical Info / Product Notes	Note: Product is not sterile.

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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)