

# CP-31398

## p53 activator

CP-31398 restores mutant p53 tumor suppressor function *in vitro* without affecting p53 homologs p63 and p73, and *in vivo* inhibiting UVB-induced skin carcinogenesis in mice. Stabilization of p53 inhibits ubiquitination without altering phosphorylation at Ser<sup>12</sup> or Ser<sup>20</sup> or MDM2 binding. Administration of this or similar compounds has the potential to enhance cytotoxic chemotherapy.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-GR360-0010	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	N'-[2-[2-(4-Methoxyphenyl)ethenyl]-4-quinazoliny]-N,N-dimethyl-1,3-propanediamine dihydrochloride
<b>Appearance</b>	Yellow solid.
<b>CAS</b>	259199-65-0
<b>Couple Target</b>	p53
<b>Couple Type</b>	Activator
<b>Formula</b>	$C_{22}H_{26}N_4O \cdot 2HCl$
<b>MW</b>	435.4
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in water (25mg/ml) or DMSO (5mg/ml).
<b>Source</b>	Synthetic.

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-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)