

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 Ser¹² or Ser²⁰ or MDM2 binding. Administration of this or similar compounds has the potential to enhance cytotoxic chemotherapy.

Citations: 5

[View Online »](#)

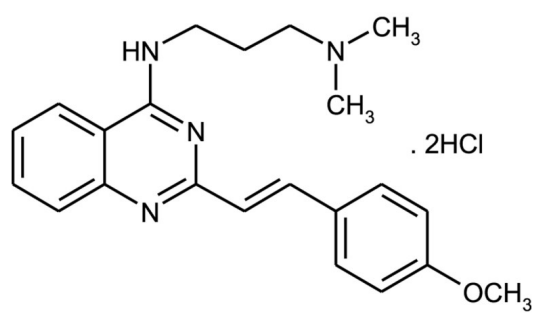
Ordering Information

[Order Online »](#)

BML-GR360-0010	10mg
BML-GR360-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

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

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)