

Pifithrin- α (cyclic) . hydrobromide

p53 inhibitor

Stable analog of pifithrin- α (Prod. No. BML-GR325), with biological properties similar to its parent compound, but with reduced cytotoxicity. Inhibitor of p53. Reversibly inhibits p53-dependent transactivation of p53-responsive genes and reversibly blocks p53-mediated apoptosis. Potent STAT6 transcriptional inhibitor. Increases apoptosis in target cells when used in combination with antimicrotubule agents.

Citations: 4

[View Online »](#)

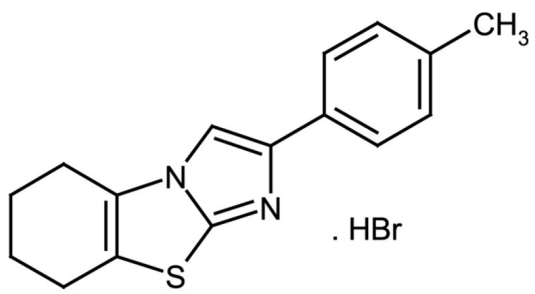
Ordering Information

[Order Online »](#)

ALX-270-287-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	60477-34-1
Couple Target	p53
Couple Type	Inhibitor
Formula	$C_{16}H_{16}N_2S \cdot HBr$
MW	268.4 . 80.9
Purity	≥95% (HPLC, TLC)
Solubility	Soluble in DMSO (100mg/ml).

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)