

Geldanamycin

HSP90 inhibitor

Potent antitumor antibiotic. Inhibitor of pp60src tyrosine kinase and of c-myc gene expression in murine lymphoblastoma cells. Inhibits the transforming activity of abl, erbB, fps, src, and yes. Binds specifically to heat shock protein 90 (HSP90) and to its endoplasmic reticulum homolog GP96 (GRP94). Capable of destabilizing several oncogene and proto-oncogene products. Potent inhibitor of the nuclear hormone receptor family. Protects against α -synuclein toxicity to dopaminergic neurons in Drosophila. Destabilizes mutant p53 protein from a number of breast, leukemic, and prostate cell lines. Inhibits basal and hypoxia-induced expression of c-Jun (IC_{50} =75nM) and abolishes hypoxia-induced increase in c-Jun N-terminal kinase (JNK) activity. Inhibits telomerase activity through inhibition of HSP90, a chaperone required for the assembly and activation of telomerase in human cells. Hsp90 inhibitors geldanamycin or 17-DMAG reduces the uptake of Chaperone Mediated Autophagy (CMA) substrates by isolated lysosomes. ~10-fold more potent than herbimycin A (Prod. No. BML-EI227)

Citations: 39

[View Online »](#)

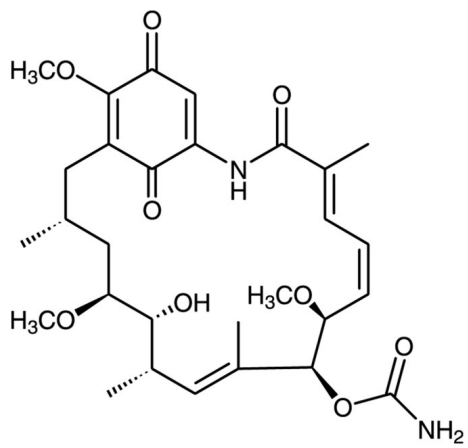
Ordering Information

[Order Online »](#)

BML-EI280-0001	1mg
----------------	-----

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.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	30562-34-6
Couple Target	HSP90, Src kinase
Couple Type	Inhibitor
Formula	$C_{29}H_{40}N_2O_9$
MW	560.6
Purity	≥95% (HPLC)
RTECS	LX8920000
Solubility	Soluble in DMSO (10mg/ml); insoluble in water.
Technical Info / Product Notes	Replacement for ADI-HPK-102

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)