

17-AAG

HSP90 inhibitor

17-AAG is a less toxic and more stable analog of geldanamycin. It is an HSP90 inhibitor that displays a 100-fold higher affinity for HSP90 derived from tumor cells compared to HSP90 from normal cells. 17-AAG inhibits Akt activation and expression in tumors and synergizes with a number of antitumor agents such as taxol, cisplatin, and UCN-014. 17-AAG causes the inactivation, destabilization and eventual degradation of HIF-1 α . Inhibitor of telomerase activity. Inducer of apoptosis with antitumor activity. Inducer of macroautophagy.

Citations: 37

[View Online »](#)

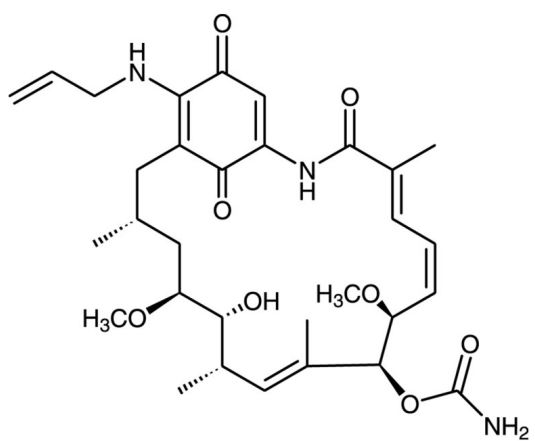
Ordering Information

[Order Online »](#)

BML-EI308-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	17-(Allylamino)-17-desmethoxygeldanamycin
Appearance	Red to dark red or purple solid.
CAS	75747-14-7
Couple Target	HSP90
Couple Type	Inhibitor
Formula	$C_{31}H_{43}N_3O_8$
Identity	Determined by ¹ H-NMR.
MW	585.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>20mg/ml) or 100% ethanol (10mg/ml).
Source	Semisynthetic derivative from geldanamycin.
Technical Info / Product Notes	Replacement for ADI-HPK-101 .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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