

# 17-AAG

17-Allylamino-17-demethoxygeldanamycin; Heat shock protein 90 (Hsp90) inhibitor

Catalog # ant-agl-5, ant-agl-25

For research use only

Version # 16D09-MM

## PRODUCT INFORMATION

### Content:

17-AAG is supplied as a lyophilized purple powder. It is available in two pack sizes:

- **ant-agl-5:** 5 mg
- **ant-agl-25:** 25 mg

### Storage and stability:

- 17-AAG is shipped at room temperature. Store lyophilized product at -20°C. Protect from light. Lyophilized product is stable for 6 months when stored properly.

- Once solubilized 17-AAG should be aliquoted and stored in the dark at -20°C for 3 months. Avoid repeated freeze-thaw cycles. Protect from light. Resuspended product is stable for 3 months when properly stored.

## QUALITY CONTROL

- Purity: ≥95% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) is confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## BACKGROUND

17-Allylamino-17-demethoxygeldanamycin (17AAG) is a less toxic and more stable analogue of geldanamycin (GA)<sup>1</sup>. Even though 17-AAG binding to heat shock protein 90 (Hsp90) is weaker than GA, 17-AAG displays similar antitumor effects as GA and a better toxicity profile.

17-AAG is currently in phase I clinical trial in several centers worldwide. Preliminary data obtained from these trials demonstrate that antitumor activity is achieved at concentrations below the maximum tolerated dose<sup>2</sup>.

**1. Schulte TW. & Neckers LM., 1998.** The benzoquinone ansamycin 17-allylamino-17-demethoxygeldanamycin binds to HSP90 and shares important biologic activities with geldanamycin. *Cancer Chemother Pharmacol* 42(4):273-9. **2. Agnew EB. et al. 2001.** Measurement of the novel antitumor agent 17-(allylamino)-17-demethoxygeldanamycin in human plasma by high-performance liquid chromatography. *J Chromatogr B Biomed Sci Appl* 755:237-43.

## CHEMICAL PROPERTIES

**CAS number:** 75747-14-7

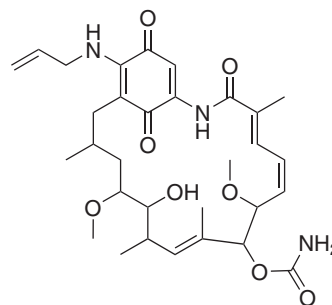
**Formula:** C<sub>31</sub>H<sub>43</sub>N<sub>3</sub>O<sub>8</sub>

**Molecular weight:** 586

**Solubility:** Soluble in DMSO (10 mg/ml)

**Working concentration:** 10 nM - 10 μM

### Structure:



## RELATED PRODUCTS

Product	Catalog Code
Geldanamycin	ant-gl-5
FITC-Geldanamycin	ant-fgl-1
17-DMAG (water soluble GA analogue)	ant-dgl-5

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Hong Kong: +852 3-622-34-80  
E-mail: info@invivogen.com