

# 17-DMAG

Water-soluble geldanamycin analogue

Catalog # ant-dgl-5, ant-dgl-25

For research use only

Version # 16G04-MM

## PRODUCT INFORMATION

### Contents:

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

- **ant-dgl-5:** 5 mg
- **ant-dgl-25:** 25 mg

### Storage and stability:

- 17-DMAG is shipped at room temperature. Store lyophilized product at -20°C. Protect from light. Lyophilized 17-DMAG is stable for 6 months when stored properly.
- Once solubilized 17-DMAG should be aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles. Protect from light. Resuspended product is stable for 3 months when properly stored.

## QUALITY CONTROL

- Purity: ≥92% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) is confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

17-(Dimethylaminoethylamino)-17-demethoxygeldanamycin (17-DMAG, also known as alvespimycin) is the first water-soluble analog of 17-AAG. This Hsp90 inhibitor shows promise in preclinical models<sup>1</sup>. 17-DMAG has excellent bioavailability, is widely distributed to tissues, and is quantitatively metabolized much less than is 17-AAG<sup>2</sup>. Clinical studies have shown that alvespimycin (KOS-1022) has therapeutic effects and is well tolerated in patients with solid tumours, advanced solid tumours or acute myeloid leukaemia, indicating that it has potential to be a new drug of choice for cancer therapy<sup>2</sup>.

**1. Workman P., 2003.** Overview: translating Hsp90 biology into Hsp90 drugs. *Curr Cancer Drug Targets* 3(5):297-300. **2. Egorin MJ, et al., 2002.** Pharmacokinetics, tissue distribution, and metabolism of 17-(dimethylaminoethylamino)-17-demethoxygeldanamycin (NSC 707545) in CD2F1 mice and Fischer 344 rats. *Cancer Chemother Pharmacol* 49(1):7-19.

## CHEMICAL PROPERTIES

**CAS number:** 467214-21-7

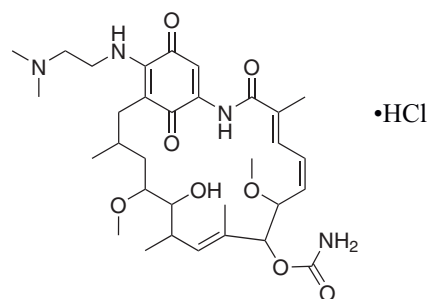
**Synonyms:** NSC 707545, Alvespimycin HCl

**Formula:** C<sub>32</sub>H<sub>48</sub>N<sub>4</sub>O<sub>8</sub> •HCl

**Molecular weight:** 653.2

**Solubility:** H<sub>2</sub>O (10 mg/ml)

**Structure:**



The use of 17-DMAG is covered under US Patent 6,890,917 owned and licensed by the NIH to InvivoGen.

## RELATED PRODUCTS

Product	Catalog Code
17-AAG (GA analogue)	ant-agl-5
FITC-Geldanamycin	ant-fgl-1
Geldanamycin	ant-gl-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