

17-DMAG

Water-soluble geldanamycin analogue

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

For research use only

Version # 09E25-MM

PRODUCT INFORMATION

Contents:

17-DMAG is supplied as a lyophilized purple powder:

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

Storage and stability:

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

Quality control

Purity controlled by TLC: 99%

DESCRIPTION

17-(Dimethylaminoethylamino)-17-demethoxygeldanamycin (17-DMAG, NSC 707545) is the first water-soluble analogue of 17-AAG. This Hsp90 inhibitor shows promise in preclinical models¹. 17-DMAG has excellent bioavailability, is widely distributed to tissues, and is quantitatively metabolized much less than is 17-AAG².

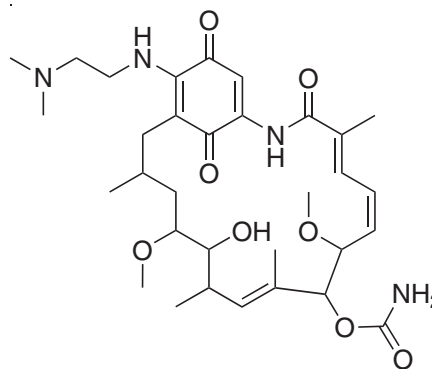
1. Workman P., 2003. Overview: translating Hsp90 biology into Hsp90 drugs. *Curr Cancer Drug Targets* 3(5):297-300. **2. Egorin MJ, et al., 2002.** Pharmacokinetics, tissue distribution, and metabolism of 17-(dimethylaminoethylamino)-17-demethoxygeldanamycin (NSC 707545) in CD2F1 mice and Fischer 344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

CHEMICAL PROPERTIES

Formula: C₃₂H₄₈N₄O₈, HCl

Molecular weight: 617

Solubility: H₂O (10 mg/ml).



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
Geldanamycin	ant-gl-5
17-AAG (GA analogue)	ant-agl-5
17-AEP-GA (water soluble GA analogue)	ant-egl-5
17-DMAP-GA (water soluble GA analogue)	ant-mgl-5
Biotin-GA	ant-bgl-1

TECHNICAL SUPPORT

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
Europe: +33 562-71-69-39
E-mail: info@invivogen.com
Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite A
San Diego, CA 92121 - USA