

# 17-DMAP-GA

Water-soluble geldanamycin analogue

Catalog # ant-mgl-1, ant-mgl-5

For research use only

Version # 09E25-MM

## PRODUCT INFORMATION

### Content:

17-DMAP-GA is supplied as a lyophilized purple powder:

-ant-mgl-1: 1 mg

-ant-mgl-5: 5 mg

### Storage and stability:

- 17-DMAP-GA is shipped at room temperature. Store lyophilized 17-DMAP-GA at -20°C in the dark. Lyophilized 17-DMAP-GA is stable for 6 months when stored properly.

- Once solubilized 17-DMAP-GA should be aliquoted and stored in the dark at -20°C for 6 months. Avoid repeated freeze-thaw cycles. Resuspended product is stable for 3 months when properly stored.

### Quality control:

Purity controlled by TLC 99%.

## BACKGROUND

17-(Dimethylaminopropylamino)-17-demethoxygeldanamycin (17-DMAP-GA) belongs to a new set of geldanamycin analogues that have been synthesized based on binding affinity to Hsp90 and water solubility. 17-DMAP-GA was shown to greatly inhibit the growth of cancer cells (IC<sub>50</sub> below 100 nM)<sup>1</sup>. Its binding affinity to Hsp90 was not significantly affected while its water solubility was highly improved compared to 17-AAG.

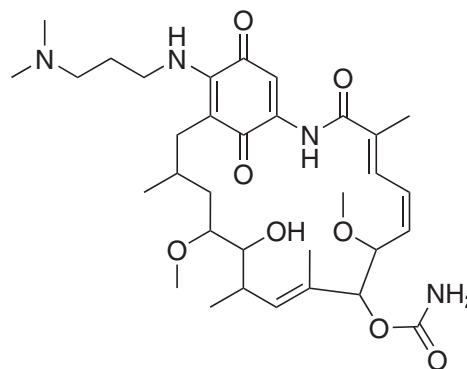
1. Tian ZQ. *et al.*, 2004. Synthesis and biological activities of novel 17-aminogeldanamycin derivatives. *Bioorg Med Chem.* 12(20):5317-29.

## CHEMICAL PROPERTIES

**Formula:** C<sub>33</sub>H<sub>50</sub>N<sub>4</sub>O<sub>8</sub>

**Molecular weight:** 631

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



## RELATED PRODUCTS

Product	Catalog Code
Geldanamycin	ant-gl-5
17-AAG (GA analogue)	ant-agl-5
17-DMAG (water soluble GA analogue)	ant-dgl-5
17-AEP-GA (water soluble GA analogue)	ant-egl-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](mailto:info@invivogen.com)

Website: [www.invivogen.com](http://www.invivogen.com)



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