

17-AEP-GA

Water-soluble geldanamycin analogue

Catalog # ant-egl-1, ant-egl-5

For research use only

Version # 09E26-MM

PRODUCT INFORMATION

Contents:

17-AEP-GA supplied as a lyophilized purple powder:

-ant-egl-1: 1 mg

-ant-egl-5: 5 mg

Storage and stability:

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

- Once solubilized 17-AEP-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-[2-(Pyrrolidin-1-yl)ethyl]amino-17-demethoxygeldanamycin (17-AEP-GA) is a new geldanamycin (GA) analogue with an alkylamino group in place of the methoxy moiety at C17. 17-AEP-GA is less cytotoxic than GA and remains biologically active¹. 17-AEP-GA was shown to induce similar tumor cell growth inhibition than 17-AAG and, unlike 17-AAG which is soluble in DMSO, to be water soluble.

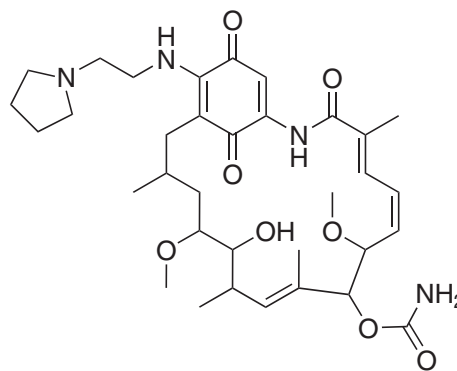
1. Tian ZQ. *et al.*, 2004. Synthesis and biological activities of novel 17-aminogeldanamycin derivatives. *Bioorg Med Chem.* 12(20):5317-29.

CHEMICAL PROPERTIES

Formula: C₃₄H₅₀N₄O₈

Molecular weight: 643

Solubility: H₂O (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-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