

Geldanamycin

An inhibitor of the heat-shock protein Hsp 90

Catalog ant-gl-5, ant-gl-25

For research use only

Version # 09E25-MM

PRODUCT INFORMATION

Content:

Geldanamycin is supplied as a yellow powder.

- **ant-gl-5:** 5 mg
- **ant-gl-25:** 25 mg

Storage and stability:

- Geldanamycin is shipped at room temperature. Store lyophilized geldanamycin at -20°C in the dark. Lyophilized geldanamycin is stable for 6 months when stored properly.
- Once solubilized geldanamycin 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%

BACKGROUND

Geldanamycin (GA) is a natural product produced by *Streptomyces hygroscopicus*. InvivoGen produces GA from a mutant strain of *S. hygroscopicus*, inactivated for the synthesis of nigericin, a common contaminant of GA. GA binds with high affinity into the ATP binding pocket of Hsp90. Hsp90 is a ubiquitous molecular chaperone critical for the folding, assembly and activity of multiple mutated and overexpressed signaling proteins that promote the growth and/or survival of tumor cells. Hsp90 client proteins include mutated p53, Raf-1, Akt, ErbB2 and hypoxia-inducible factor 1a (HIF-1a)¹. Binding of GA to Hsp90 causes the destabilization and degradation of its client proteins².

1. Neckers L., 2002. Hsp90 inhibitors as novel cancer chemotherapeutic agents. Trends Mol Med 8(4 Suppl):S55-61. 2. Whitesell L. et al., 1994. Inhibition of heat shock protein HSP90-pp60v-src heteroprotein complex formation by benzoquinone ansamycins: essential role for stress proteins in oncogenic transformation. PNAS 91(18):8324-8.

CHEMICAL PROPERTIES

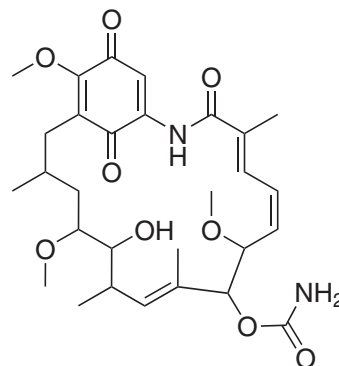
CAS n°: 30562-34-6

Formula: C₂₉H₄₀N₂O₉

Molecular weight: 560

Solubility: Soluble in DMSO (1 mg/ml)

Working concentration: 1 nM - 10 μM



RELATED PRODUCTS

Product	Catalog Code
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
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