

## Geldanamycin

Geldanamycin is an antibiotic and potent antagonist of the heat shock protein 90 (Hsp90) protein (KD=1.2uM), binding its ATP site and inhibiting its chaperone activity. Induces degradation of mutated proteins v-Src, Bcr-Abl, Raf, p53, and steroid receptors in tumor cells selectively over non-mutated counterparts via Hsp90 inhibition. Affects multiple cellular processes, including gene expression, cell proliferation, apoptosis, and angiogenesis.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1235</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 mg per vial

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	560.64 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 98%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	10 mM
<b>Reconstitution Volume:</b>	0.178 mL
<b>Reconstitution Solvent:</b>	DMSO

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)