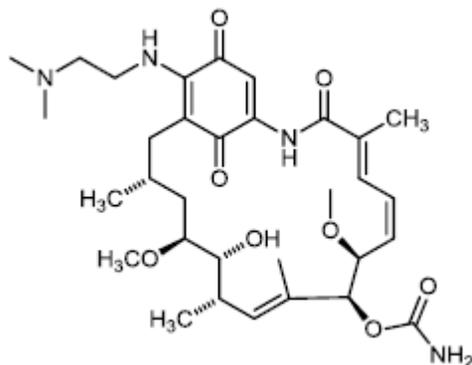


## 17-DMAG. HCl

Cat. # AC-2006

**Background:** Alvespimycin (17-DMAG) is a potent HSP90 inhibitor with  $IC_{50}$  of 62 nM in a cell-free assay. Geldanamycin analog that displays superior pharmacological properties and induces apoptosis in a variety of tumor cell lines. Inhibits angiogenesis. Cell permeable.

**Application:** Autophagy inducer



### Product Information

|                        |  |
|------------------------|--|
| <b>Purity:</b>         | >98% by TLC  |
| <b>MW:</b>             | 653.2  |
| <b>Formula:</b>        | $C_{32}H_{48}N_4O_8 \cdot HCl$   |
| <b>Physical State:</b> | Lyophilized purple powder  |
| <b>Amount:</b>         | 1mg or 5mg   |
| <b>Solubility:</b>     | 30 mg/ml in DMSO; 10 mg/mL in ethanol  |
| <b>Storage:</b>        | Store as supplied at $-20^{\circ}C$ for up to 2 years. Store solutions at $-20^{\circ}C$ for up to 3 months. |

### References

1. M Hollingshead *et al.* *Cancer Chemother. Pharmacol.* 2005 56:637
2. G Kaur *et al.* *Clin. Cancer Res.* 2004 10:4813

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2013 LifeSensors, Inc. All Rights Reserved