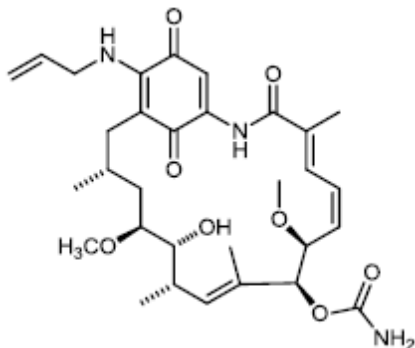


17-AAG

Cat. # AC-2005

Background: Tanespimycin (17-AAG) is a semi-synthetic analog of geldanamycin which is less toxic and more stable¹. Selectively binds to and inhibits HSP90 from tumor cells displaying a 100-fold higher affinity for tumor cell-derived over normal cell HSP90². Displays anti-angiogenic activity³. Sensitizes tumor cells to apoptosis in response to such agents as taxol⁴ and cisplatin⁵. Cell permeable.

Application: Autophagy inducer



Product Information

Purity:	>98% by TLC
MW:	585.7
Formula:	C ₃₁ H ₄₃ N ₃ O ₈
Physical State:	Lyophilized red to dark red powder
Amount:	1mg or 5mg
Solubility:	>50 mg/ml in DMSO; 5 mg/mL in ethanol
Storage:	Store desiccated at -20°C. Store solutions at -20°C for 3 months.

References

1. TW Schulte *et al.* *Cancer Chemother. Pharmacol.* 1998 42:273
2. A Kamal *et al.* *Nature* 2003 425:407
3. G Kaur *et al.* *Clin. Cancer Res.* 2004 10:4813
4. DB Solit *et al.* *Cancer Res.* 2003 63:2139
5. IA Vasilevskaya *et al.* *Mol. Pharmacol.* 2004 65:235

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