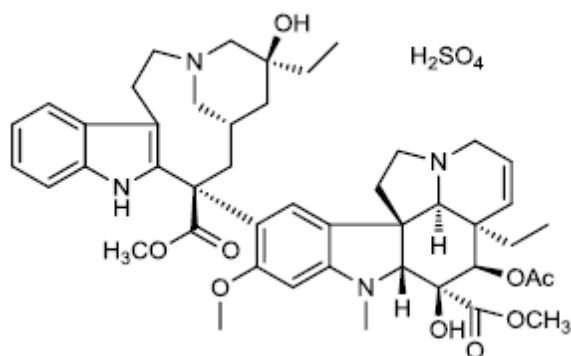


## Vinblastine sulfate

Cat. # AC-1007

**Background:** Semisynthetic alkaloidal anticancer agent. Induces cell cycle arrest at G2/M phase by inhibiting mitotic spindle formation<sup>1</sup>. Inhibits normal microtubule assembly and induces aberrant tubulin polymerization<sup>1</sup>. Induces apoptosis in a variety of cell lines<sup>2</sup>. Inhibits autophagosome maturation<sup>3</sup>. Inhibits palmitoylation of tubulin<sup>4</sup>. Vinblastine sulfate inhibits microtubule formation and suppresses nAChR activity with IC<sub>50</sub> of 8.9 μM in a cell-free assay, used to treat certain kinds of cancer.

**Application:** Autophagy inhibitor: inhibits autophagosome maturation



### Product Information

<b>Purity:</b>	98% by TLC
<b>MW:</b>	909.1
<b>Formula:</b>	C <sub>46</sub> H <sub>58</sub> N <sub>4</sub> O <sub>9</sub> · H <sub>2</sub> SO <sub>4</sub>
<b>CAS No.</b>	143-67-9
<b>Physical State:</b>	Lyophilized white crystalline powder
<b>Amount:</b>	5mg, 25mg
<b>Solubility:</b>	25 mg/mL in DMSO; 50 mg/mL in water
<b>Storage:</b>	Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month.

### References

1. E Hamel *Med. Res. Rev.* 1996 16:207
2. K Tsukidate *et al. Am. J. Pathol.* 1993 143:918
3. CP Satori *et al. Anal. Chem.* 2013 85:11391
4. JM Caron and M Herwood *Chemotherapy* 2007 53:51

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