

## NMS-873

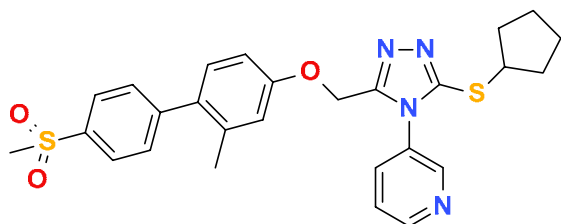
Cat. # SI9752

**Background:** p97 (aka VCP, Cdc48) is intimately involved in the ERAD process through the retrotranslocation of misfolded proteins targeted for degradation by the proteasome. It is a homohexamer of AAA<sup>+</sup> ATPase subunits and ATP hydrolysis is required for protein extraction from the ER. NMS-873 is an allosteric inhibitor of ATPase activity, binding to the D1 and D2 domains of adjacent subunits (Magnaghi, D'Alessio et al. 2013). NMS-873 is reported to have an IC<sub>50</sub> for p97 of 20-30nM.

**Application:** Inhibition of ERAD, induction of apoptosis in cancer cells

### Product Information

<b>CAS No.:</b>	1418013-75-8
<b>Purity:</b>	99%, NMR is consistent with structure
<b>Molecular Weight:</b>	520.67
<b>Physical State:</b>	Powder
<b>Quantity:</b>	5mg
<b>Solubility:</b>	192mM in DMSO
<b>Storage:</b>	Store dessicated as supplied at -20°C for 2 years
<b>Formula:</b>	C <sub>27</sub> H <sub>28</sub> N <sub>4</sub> O <sub>3</sub> S <sub>2</sub>
<b>SMILES String:</b>	<chem>Cc1cc(ccc1c2ccc(cc2)S(=O)(=O)C)OCc3nnc(n3c4cccnc4)SC5CCCC5</chem>



### References

Magnaghi, P., R. D'Alessio, et al. (2013). "Covalent and allosteric inhibitors of the ATPase VCP/p97 induce cancer cell death." *Nat Chem Biol* **9**(9): 548-556.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2015 LifeSensors, Inc. All Rights Reserved