

## NEDD8-activating enzyme inhibitor

1-[(1S,2S,4R)-4-[4-[2,3-dihydro-1H-inden-1S-ylamino]-7H-pyrrolo[2,3-d]pyrimidin-7-yl]-2-hydroxycyclopentyl] methyl sulfamate

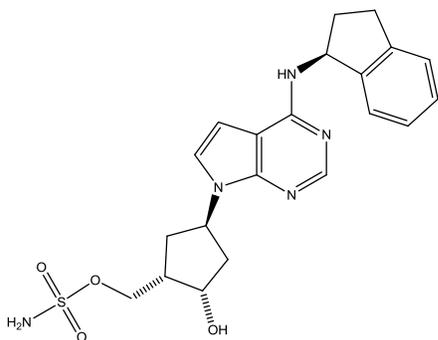
Cat. # SI9830

**Background:** Contains the active ingredient of MLN4924, a substrate-assisted inhibitor of the NEDD8-activating enzyme (NAE). NAE activates Nedd8 (Neural precursor cell expressed, developmentally down-regulated 8) for conjugation to a variety of ubiquitin E3 ligases. MLN4924 selectively inhibits NAE activity and represents a new approach to targeting the UPS for the treatment of cancer. Treatment of cells with MLN4924 results in inhibition of the NEDD8 pathway. Cell permeable.

**Application:** Inhibition of NEDD8-dependent pathways.

### Product Information

<b>CAS No.</b>	905579-51-3
<b>Purity:</b>	> 98% (TLC); NMR (Conforms)
<b>Molecular Weight:</b>	443.2
<b>Physical State:</b>	White powder
<b>Quantity:</b>	1mg
<b>Solubility:</b>	DMSO (10 mg/ml)
<b>Storage:</b>	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 3 months.



**Formula:** C<sub>21</sub>H<sub>25</sub>N<sub>5</sub>O<sub>4</sub>S

### References

1. Brownell JE et al. Substrate-assisted inhibition of ubiquitin-like protein-activating enzymes: the NEDD8 E1 inhibitor MLN4924 forms a NEDD8-AMP mimetic *in situ*. *Molecular Cell* 2010, **37**, 102-111.
2. Milhollen MA et al. MLN4924, a NEDD8-activating enzyme inhibitor, is active in diffuse large B-cell lymphoma models: rationale for treatment of NF-κB-dependent lymphoma. *Blood*, 2010, **116**,1515-1523.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2011 LifeSensors, Inc. All Rights Reserved