

## SMER3

### 9H-Indeno[1,2-e][1,2,5]oxadiazolo[3,4-b]pyrazin-9-one

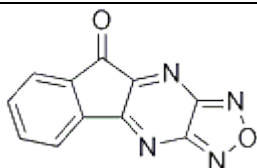
Cat. # SI9880

**Background:** Small Molecule Enhancer of Rapamycin 3 (SMER3) is a specific inhibitor of the ubiquitin E3 ligase SCF<sup>MET30</sup>. SMER3 binds directly to the F-box subunit Met30 and blocks binding to the SCF core complex. This inhibitor is cell permeable. Significant inhibition is typically seen at 30-100µM.

**Application:**

**Product Information:**

<b>CAS No.</b>	67200-34-4
<b>Purity:</b>	>98% (TLC); NMR (Conforms)
<b>Molecular Weight:</b>	224.2 Da
<b>Physical State:</b>	Yellow Powder
<b>Quantity:</b>	10mg, 50mg
<b>Solubility:</b>	May be dissolved in DMSO (20mg/ml, warm)
<b>Storage:</b>	Store desiccated as supplied at ambient temp. for up to 2 years. Store solutions at -20°C for up to 2 months.



**Formula:**

C<sub>11</sub>H<sub>4</sub>N<sub>4</sub>O<sub>2</sub>S

**References**

1. Aghajany, M. *et al.* Chemical genetics of TOR identifies an SCF family E3 ubiquitin ligase inhibitor. *Nat. Biotech.* 2010, 28(7):738.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2014 LifeSensors, Inc. All Rights Reserved