

## Epoxomicin/BU-4061T

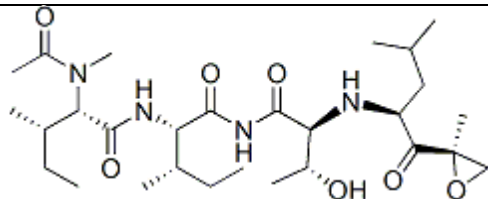
Cat. # SI9780

**Background:** One of the most potent and selective proteasome inhibitors, Epoxomicin inhibits all three activities of the proteasome. Epoxomicin (100 nM) results in a 30-fold increase in the levels of p53 protein, a known target of the proteasome, in HUVECs. Epoxomicin (10  $\mu$ M) results in the accumulation of ubiquitinated proteins in HeLa cells. Epoxomicin (10  $\mu$ M) inhibits I $\kappa$ B $\alpha$  degradation by 10-fold in HeLa cells<sup>1</sup>.

**Application:** Cell-based and in vitro studies of proteasome function

### Product Information:

<b>CAS No.</b>	134381-21-8
<b>Purity:</b>	>99.0% by HPLC
<b>Molecular Weight:</b>	554.72 Da
<b>Physical State:</b>	Powder
<b>Quantity:</b>	100 $\mu$ g
<b>Solubility:</b>	DMSO (10 mg/mL)
<b>Storage:</b>	Store desiccated as supplied at -20°C for 2 years.



**Formula:** C<sub>28</sub>H<sub>50</sub>N<sub>4</sub>O<sub>7</sub>

### References

1. Meng, L., R. Mohan, et al. (1999). "Epoxomicin, a potent and selective proteasome inhibitor, exhibits in vivo antiinflammatory activity." *Proc Natl Acad Sci U S A* **96**(18): 10403-10408.
2. Hakim, V., Cohen, L.D., et al. (2016) "The effects of proteasomal inhibition on synaptic proteostasis." *EMBO J* **35**(20): 2238-2262.

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