

## PS-341 (Pyz-Phe-boroLeu; MG-341)

### N-(2-pyrazine)carbonyl-L-phenylalanyl-L-leucyl boronic acid

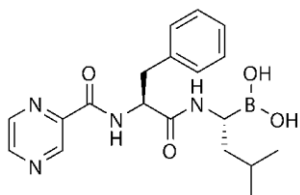
Cat. # SI9720

**Background:** The active ingredient in Bortezomib, Pyz-Phe-boroLeu is a dipeptide boronic acid inhibitor of proteasome activity ( $K_i=0.6$  nM)

**Application:** Inhibition of 26S proteasome chymotrypsin-like activity.

#### Product Information:

<b>CAS No.</b>	179324-69-7
<b>Purity:</b>	>98% (HPLC)
<b>Molecular Weight:</b>	384.24 Da
<b>Physical State:</b>	White Powder
<b>Quantity:</b>	5mg, 25mg
<b>Solubility:</b>	DMSO
<b>Storage:</b>	Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to one week.



**Formula:**



#### References

1. Yi-He Ling et al. PS-341, a Novel Proteasome Inhibitor, Induces Bcl-2 Phosphorylation and Cleavage in Association with G2-M Phase Arrest and Apoptosis. *Mol Cancer Ther* 2002;1:841-849.
2. Adams J. Development of the Proteasome Inhibitor PS-341. *The Oncologist* 2002;7:9-16.
3. Dudek SA et al. The clinically approved proteasome inhibitor PS-341 efficiently blocks influenza A virus and vesicular stomatitis virus propagation by establishing an antiviral state. *J Virol*. 2010 Sep;84:9439-51.

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