

## Carfilzomib/PR-171

Cat. # SI9790

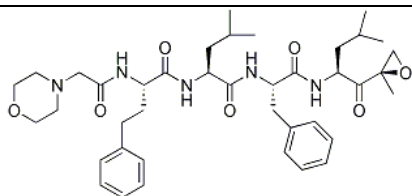
**Background:** The active ingredient in Kyprolis, carfilzomib is an epoxomicin analog proteasome inhibitor with with an IC<sub>50</sub> less than 5 nM.

**Application:**

---

**Product Information:**

<b>CAS No.</b>	868540-17-4
<b>Purity:</b>	
<b>Molecular Weight:</b>	719.91 Da
<b>Physical State:</b>	Powder
<b>Quantity:</b>	5mg
<b>Solubility:</b>	DMSO (50 mg/ml); Ethanol (<1mg/ml); Water (<1mg/ml)
<b>Storage:</b>	Store desiccated as supplied at -20°C for 2 years.



**Formula:** C<sub>40</sub>H<sub>57</sub>N<sub>5</sub>O<sub>7</sub>

---

**References**

1.

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