

# MLN2238

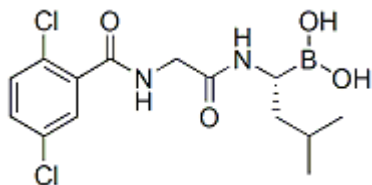
Cat. # SI9725

**Background:** MLN2238 is a proteasome inhibitor also known as ixazomib. MLN2238 is the biologically active form of MLN9708.<sup>1</sup> Under physiological conditions, the MLN9708 boronic citrate ester is rapidly hydrolyzed to a boronic acid. Much like Bortezomib, MLN2238 reversibly inhibits the trypsin- and chymotrypsin-like activities of the proteasome.<sup>2</sup>

**Application:** *In vitro* and cellular studies of proteasome function.

## Product Information:

<b>CAS No.</b>	1072833-77-2
<b>Purity:</b>	>99.0% by HPLC
<b>Molecular Weight:</b>	361.03 Da
<b>Physical State:</b>	Powder
<b>Quantity:</b>	5 mg
<b>Solubility:</b>	DMSO (72 mg/mL); ethanol (9 mg/mL); Water (<1 mg/mL)
<b>Storage:</b>	Store desiccated as supplied at -20°C for 2 years.



**Formula:** C<sub>14</sub>H<sub>19</sub>BCl<sub>2</sub>N<sub>2</sub>O<sub>4</sub>

## References

1. Kupperman, E., Lee, E.C. et al. (2010). "Evaluation of the proteasome inhibitor MLN9708 in preclinical models of human cancer." *Cancer Res* **70**(5): 1970-1980.
2. Schrader, J., Hennenberg, F., et al. (2016). "The inhibition mechanism of human 20S proteasomes enables next-generation inhibitor design." *Science* **353**(6299): 594-598.

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