

Oprozomib/ONX0912

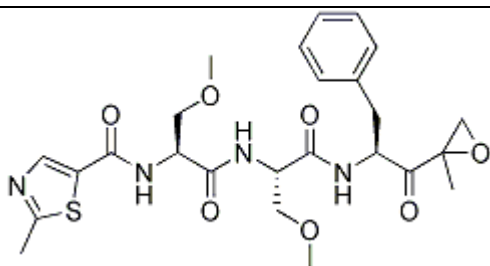
Cat. # SI9735

Background: Clinical-stage proteasome inhibitor that selectively inhibits the chymotrypsin-like activity of the 20S proteasome β 5/LMP7 with IC_{50} of 36 nM/82 nM.

Application:

Product Information:

CAS No.	935888-69-0
Purity:	>99.0% by HPLC
Molecular Weight:	532.61 Da
Physical State:	Powder
Quantity:	5mg
Solubility:	DMSO (107mg/ml); Ethanol (<1mg/ml); Water (<1mg/ml)
Storage:	Store desiccated as supplied at -20°C for 2 years.



Formula: C₂₅H₃₂N₄O₇S

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