

Product Information

Description:	(R)-MG132 is a potent, cell permeable and reversible inhibitor of proteasomes. It can inhibit proteasome activity in lysates of J558L multiple myeloma cells and EMT6 breast cancer cells. In comparison to (S)-MG132, the (R)-MG132 stereoisomer is a more effective inhibitor of chymotrypsin-like (ChTL), trypsin-like (TL), and peptidylglutamyl peptide hydrolyzing proteasome (PGPH) activities.
Chemical Formula:	C ₂₆ H ₄₁ N ₃ O ₅
Appearance:	A crystalline solid
Stability:	Up to 2 years at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
Quality Control:	The purity was determined by HPLC analysis.
CAS Registry #:	1211877-36-9
MW:	475.6 Da
Purity:	≥97%
Biological Activity:	(R)-MG132 inhibits ChTL, TL and PGPH with IC ₅₀ values of 0.22 μM, 34.4 μM, and 2.95 μM, respectively.
Solubility:	Soluble in ethanol, DMSO and DMF, purged with an inert gas. Solubility in these solvents is approximately 25 mg/ml. Do not store aqueous solutions for more than one day.

Structure

