

## MG-132 (Z-Leu-Leu-Leu-CHO)

### N-(benzyloxycarbonyl)leucinyl-leucinyl-leucinal

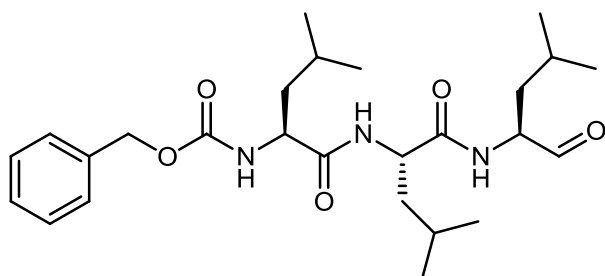
Cat. # SI9710

**Background:** MG132 is a potent, cell-permeable inhibitor of the proteasome ( $IC_{50} = 0.1 \mu\text{M}$ ) and calpain ( $IC_{50} = 1.2 \mu\text{M}$ ). MG132 inhibits TNF- $\alpha$ -induced NF- $\kappa\text{B}$  activation and I $\kappa\text{B}\alpha$  degradation ( $IC_{50} = 3 \mu\text{M}$ ). It also reduces the degradation of ubiquitin-conjugated proteins in different cell types and promotes apoptosis via activation of c-Jun N-terminal kinase (JNK1).

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

#### Product Information

<b>CAS No.</b>	133407-82-6
<b>Purity:</b>	> 98% (HPLC)
<b>Molecular Weight:</b>	475.6
<b>Physical State:</b>	White powder
<b>Quantity:</b>	10 mg, 25 mg
<b>Solubility:</b>	DMSO (45mg/ml) or Ethanol (45mg/ml)
<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:**  $\text{C}_{26}\text{H}_{41}\text{N}_3\text{O}_5$

#### References

1. Tsubuki S et al. Differential inhibition of calpain and proteasome activities by peptidyl aldehydes of di-leucine and tri-leucine. *J. Biochem.* 1996, **119**:572
2. Wu WK et al. Macroautophagy and ERK phosphorylation counteract the antiproliferative effect of proteasome inhibitor in gastric cancer cells. *Autophagy*, 2010, **6**:22.
3. Meriin AB et al. Proteasome inhibitors activate stress kinases and induce Hsp72. Diverse effects on apoptosis. *J. Biol. Chem.* 1998, **273**:6373

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