

Suc-Leu-Leu-Val-Tyr-AMC (Suc-LLVY-AMC), Proteasome Substrate

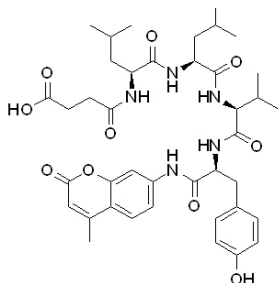
Cat. # PS500

Background: Fluorogenic substrate for measuring the chymotrypsin-like peptidase activity of the 20S proteasome, calpains and other chymotrypsin-like proteases (Ex.: 380nm Em.: 460nm).

Product Information

Quantity:	5mg
Molecular weight:	763.9
Formula:	$C_{40}H_{53}N_5O_{10}$
State:	Lyophilized powder
Purity:	>98% by TLC
Solubility:	DMSO at least 5mM. Concentration range 10-100 μ M.
Storage:	Dry reagent at 4°C; Store DMSO stock at -20°C. Avoid repeated freeze/thaw cycles

Formula:



References

1. Tsubuki S et al. (1993) Purification and characterization of a Z-Leu-Leu-Leu-MCA degrading protease expected to regulate neurite formation: a novel catalytic activity in proteasome. *Biochem. Biophys. Res. Commun.* 196, 1195
2. Stein RL. et al. (1996) Kinetic characterization of the chymotryptic activity of the 20S proteasome. *Biochem.* 35:3899-3908
3. Dang LC. et al. (1998) Kinetic and mechanistic studies on the hydrolysis of ubiquitin C-terminal 7-amido-4-methylcoumarin by deubiquitinating enzymes *Biochemistry* 37, 1868
4. Wang KK et al. (1996) An alpha-mercaptoacrylic acid derivative is a selective nonpeptide cell-permeable calpain inhibitor and is neuroprotective. *Proc. Natl. Acad. Sci. USA* 93, 6687

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved