

Background

Ubiquitin-rhodamine 110 is a quenched, fluorescent substrate for deubiquitylases, especially ubiquitin C-terminal hydrolases. Cleavage of the amide bond between the C-terminal glycine of ubiquitin and rhodamine results in an increase in rhodamine fluorescence at 535 nm (Exc. 485 nm).

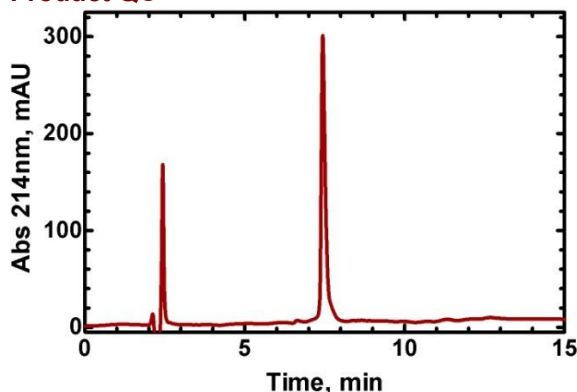
Application(s)

Measurement of deubiquitylase activity.

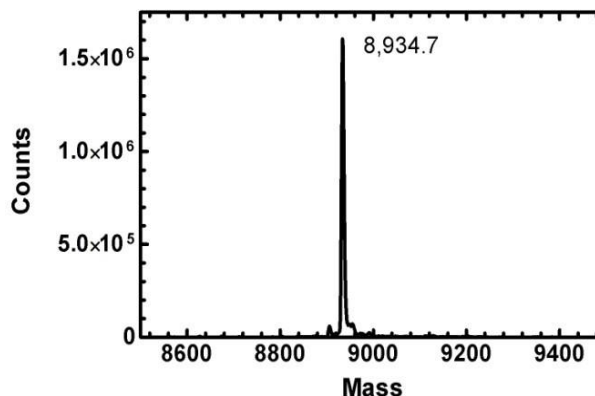
Product Specifications

Tag	None
Purity	≥ 95% by RP-HPLC
Molecular Weight	8934.2 Da
Quantity	50, 250, 500 µg
Species	Human
Expression System	<i>E. Coli</i>
Physical State	Lyophilized powder
Solubility	Aqueous buffers, DMSO
Storage	-80°C. Avoid repeated freeze/thaw cycles

Product QC



RP-HPLC



Deconvoluted MS

References

1. Tirrell DA, et al. J Biol Chem. (2013);288(25):18304-18312.
2. Greene, W. et al. (2012) PLoS Pathog. 8:e1002703.
3. Tong, X. et al. (2012) J. Biol. Chem. 287:25280.
4. Wei, W. et al. (2004) Nature 428:194.
5. Wertz, I.E. et al. (2004) Nature 430:694.
6. Hassiepen, U. et al. (2007) Anal. Biochem. 371: 201

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved