

Ubiquitin-rhodamine 110

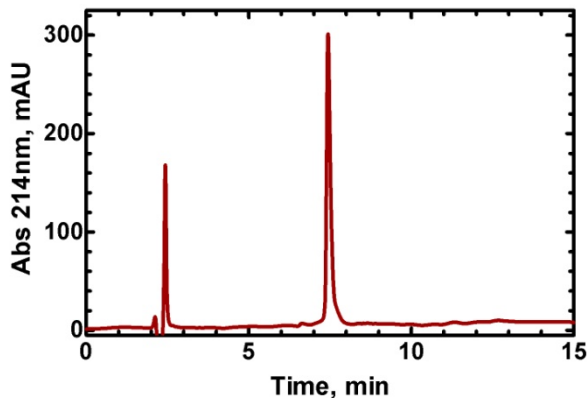
Cat. # SI230

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).

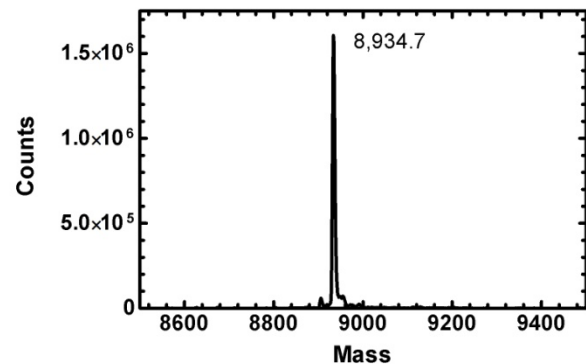
Application: Measurement of deubiquitylase activity.

Product Information

Purity:	≥ 95% by RP-HPLC
Molecular Weight:	8934.2 Da
Physical State:	Lyophilized
Quantity:	50, 250, 500 µg
Solubility:	Aqueous buffers, DMSO
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted mass spectrum

References.

- Tirat, A., Schilb, A., Riou, V., Leder, L., Gerhartz, B., Zimmermann, J., Worpenberg, S., Eidhoff, U., Freuler, F., Stettler, T., Mayr, L., Ottl, J., Leuenberger, B., and Filipuzzi, I. 2005. Synthesis and characterization of fluorescent ubiquitin derivatives as highly sensitive substrates for the deubiquitinating enzymes UCH-L3 and USP-2. *Anal. Biochem.* **343**,244-255.
- Hassiepin, U., Eidhoff, U., Meder, G., Bulber, J.F., Hein, A., Bodendorf, U., Lothois, E., and Martoglio, B., 2007. A sensitive fluorescence intensity assay for deubiquitinating proteases using ubiquitin-rhodamine 110-glycine as substrate. *Anal. Biochem.* **371**,201-207.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved