

## SUMO2 Rhodamine 110

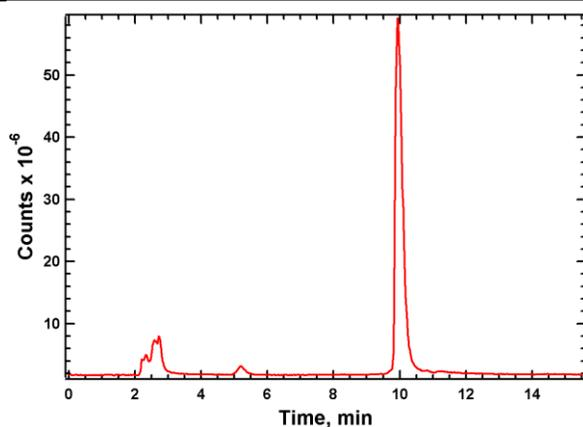
Cat. # SI530

**Background:** Human SUMO2-rhodamine 110 is a quenched, fluorescent substrate for desumoylases, especially SenP1 and SenP2. Cleavage of the amide bond between the C-terminal glycine of SUMO2 and rhodamine results in an increase in rhodamine fluorescence at 535 nm (Exc. 485 nm)..

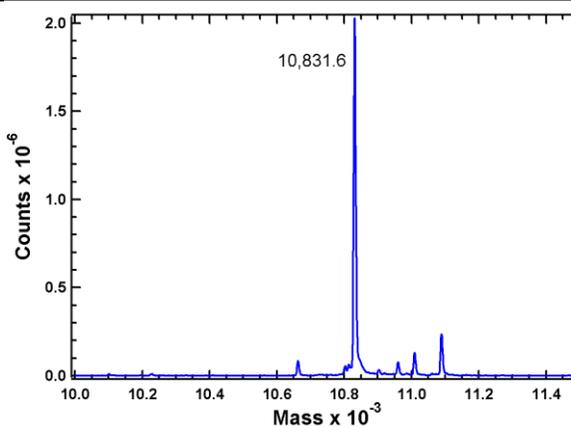
**Application:** Measurement of desumoylase activity.

### Product Information

|                          |   |
|--------------------------|---|
| <b>Purity:</b>           | ≥ 95% by RP-HPLC                          |
| <b>Molecular Weight:</b> | 10,831.1 Da                               |
| <b>Physical State:</b>   | Lyophilized powder                        |
| <b>Quantity:</b>         | 50 µg                                     |
| <b>Storage:</b>          | -80° C. Avoid repeated freeze/thaw cycles |



**RP-HPLC**



**Deconvoluted mass spectrum**

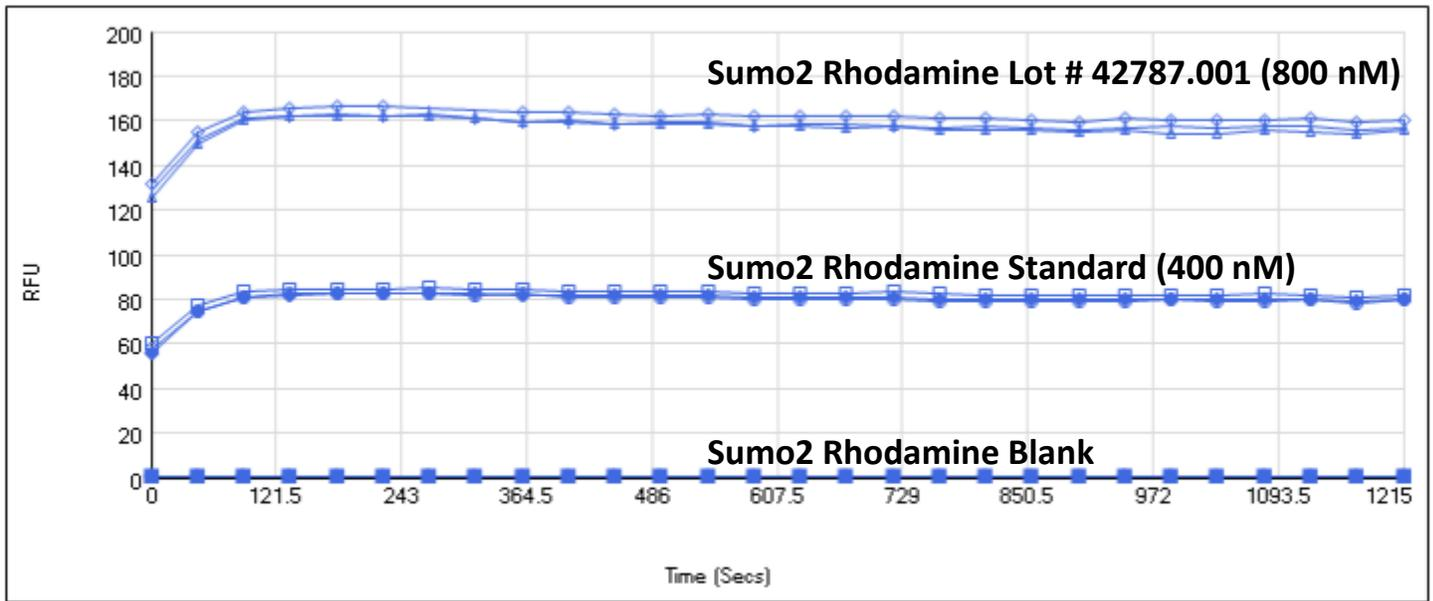
### References

1. Shanmugham A and Ovaa H. (2008) DUBs and disease: activity assays for inhibitor development. *Curr Opin Drug Discov Devel.* **11**,688-96.
2. Mason, D.E., Ek, J., Peters, E.C. and Harris J.L. (2004) Substrate profiling of deubiquitin hydrolases with a positional scanning library and mass spectrometry. *Biochemistry* **43**,6535-44.
3. Dang, L. C., F. D. Melandri, and R. L. Stein. (1998) Kinetic and mechanistic studies on the hydrolysis of ubiquitin C-terminal 7-amido-4-methylcoumarin by deubiquitinating enzymes. *Biochemistry* **37**,1868-1879.
4. Stein RL, Chen Z, Melandri F. (1995) Kinetic studies of isopeptidase T: modulation of peptidase activity by ubiquitin. *Biochemistry* **34**,12616-23.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2010 LifeSensors, Inc. All Rights Reserved

SUMO2 Rhodamine 110  
 Cat. # SI530, Lot # 42787.001

**Specific activity test in the presence of SUMO Protease**



-Lm1

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
 Copyright © 2010 LifeSensors, Inc. All Rights Reserved