

## hSUMO2-AMC

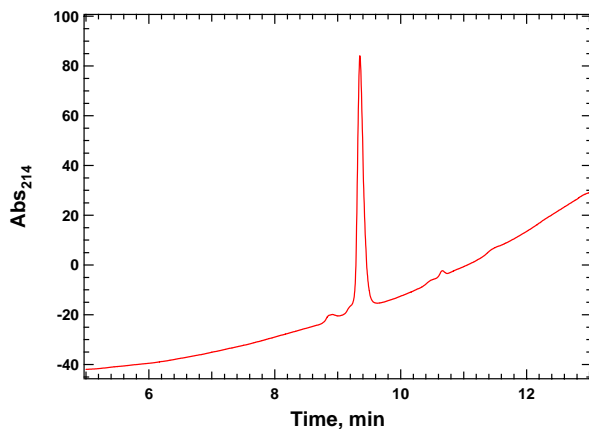
Cat. # SI520

**Background:** hSUMO2-AMC is prepared via the conjugation of 7-amino-4-methylcoumarin (AMC) to the C-terminal Gly of hSUMO2. This conjugation quenches the intrinsic fluorescence of AMC. Upon incubation with a protease recognizing hSUMOs, such as Senp1 or Senp2, AMC is released and the increase in fluorescence at 460 nm (Exc. At 380 nm) can be measured. The enhanced purity of hSUMO2-AMC compared to that of other commercial vendors leads to a greater signal to noise ratio. This protein contains no extraneous tags.

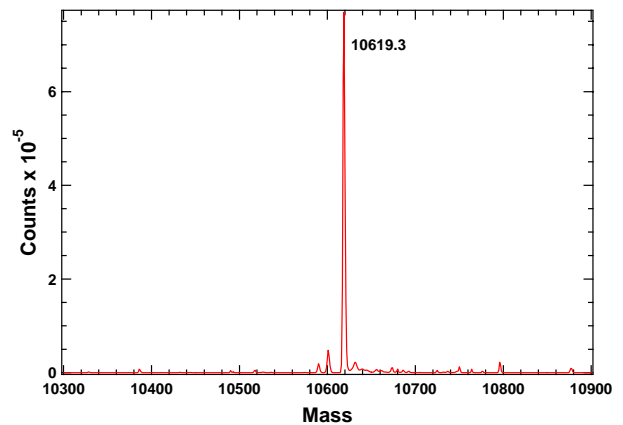
**Application:** Used for determining the activity and specificity of SUMO-specific hydrolases. High throughput screening for modulators of desumoylase activity.

### Product Information

<b>Purity:</b>	≥ 95% by RP-HPLC
<b>Molecular Weight:</b>	10,619.3 Da by MS (calculated 10,618.8 Da)
<b>Physical State:</b>	lyophilized
<b>Quantity:</b>	50 µg
<b>Solubility:</b>	
<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 © 2009 LifeSensors, Inc. All Rights Reserved