

Ubiquitin vinylmethyl ester (Ub-VME)

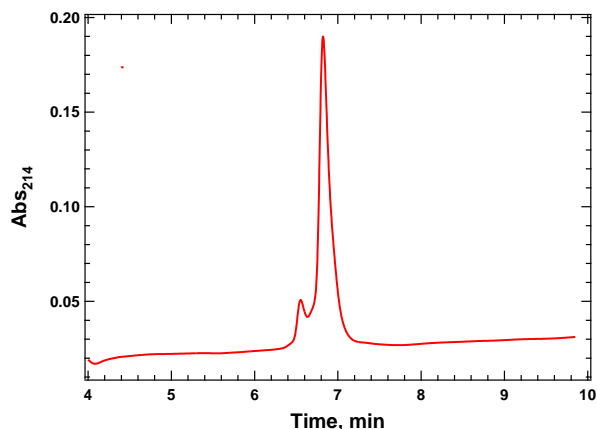
Cat. # SI240

Background: Ub-VME is synthesized by the conjugation of 4-amino-but-2-enoic methyl ester to the C-terminus of Ubiquitin Δ G76. Binding of ubiquitin by deubiquitylases (DUBs), for instance UCH's and most USPs, positions the reactive vinyl bond next to the sulfhydryl-group of the active site cysteine. Nucleophilic attack by the sulfhydryl on the vinyl bond produces a stable, covalent thioether bond between Ubiquitin and the deubiquitylase. Thus Ub-VME is a potent suicide inhibitor of DUBs.

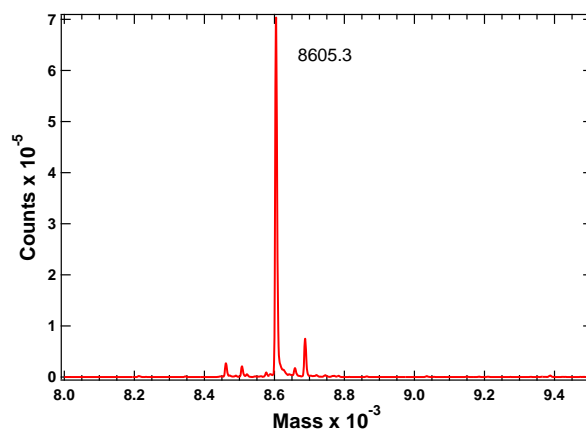
Application: This inhibitor is useful for labeling DUBs *in situ* as well as preserving the integrity of polyubiquitin chains on modified proteins for analysis or purification.

Product Information

Purity:	$\geq 95\%$ by RP-HPLC
Molecular Weight:	8,604.9 Da
Physical State:	lyophilized
Quantity:	50 μ g
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted mass spectrum

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