

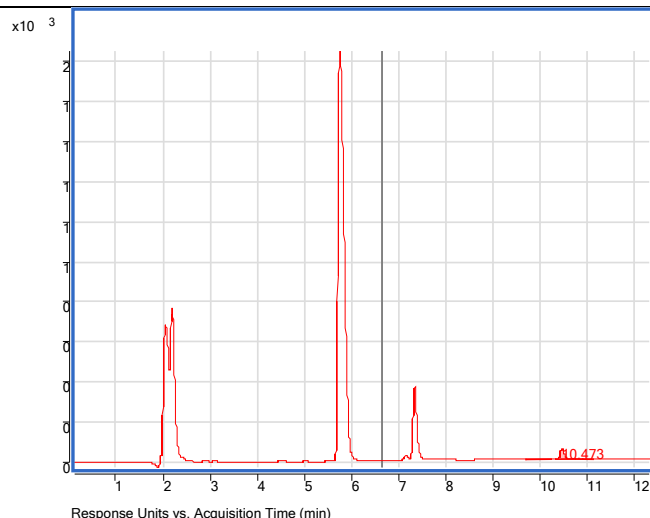
Monoubiquitinated Histone H2B Peptide (Ac-H2B[110-125](Ub)-CONH₂) Cat. # UP110

Background: Monoubiquitination of lysine 120 on human histone H2B is associated with transcriptional activity. The RNF20/40 complex functions as the E3 ligase and Ubch6 as the E2 conjugating enzyme to monoubiquitinate this site to stimulate HOX gene expression (1). Monoubiquitination of lysine 120 increases histone chaperone FACT function resulting in transcript elongation and the generation of longer transcripts (2).

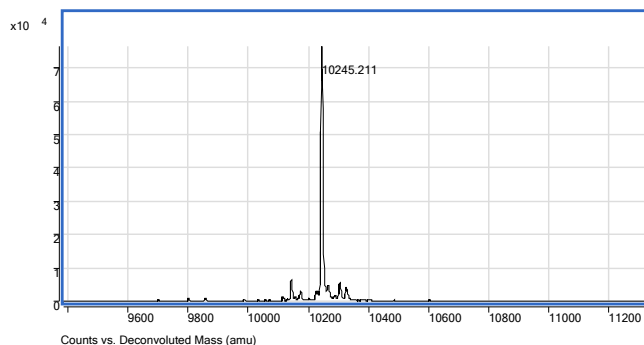
Application: This monoubiquitinated peptide can be used for deubiquitination assays or to study ubiquitin binding domain recognition.

Product Information

Sequence:	Ac-AVSEGTKAVTK(ubiquitin)YTSSK-NH ₂
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	10,244.7 Da
Physical State:	Liquid, 1mg/mL PBS
Quantity:	50µg
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted MS-spectrum

References

1. Zhu, B., et al., Monoubiquitination of human histone H2B: the factors involved and their roles in HOX gene regulation. *Mol Cell*, 2005. **20**(4):601-11.
2. Pavri, R., et al., Histone H2B monoubiquitination functions cooperatively with FACT to regulate elongation by RNA polymerase II. *Cell*, 2006. **125**(4):703-17.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved