

# Histone H2A.ZK4, K8, K12ac Peptide, Biotinylated

Catalog No. 12-0201

Lot No. 17277001

Quantity 50 µg



EpiCypher™

## Product Description:

A synthetic peptide derived from human histone H2A.Z amino acids 1-19, acetylated at lysines 4, 8, and 12. The alpha-amino group on the N-terminus is acetylated, with a PEG spacer containing a biotin moiety at the C-terminus.

## Formulation:

50 µg of lyophilized powder.

## Sequence:

Ac-AGGK(Ac)AGK(Ac)DSGK(Ac)AKTKAVSR-Peg-Biot

## Molecular Weight:

2413.8 Da

## Purity:

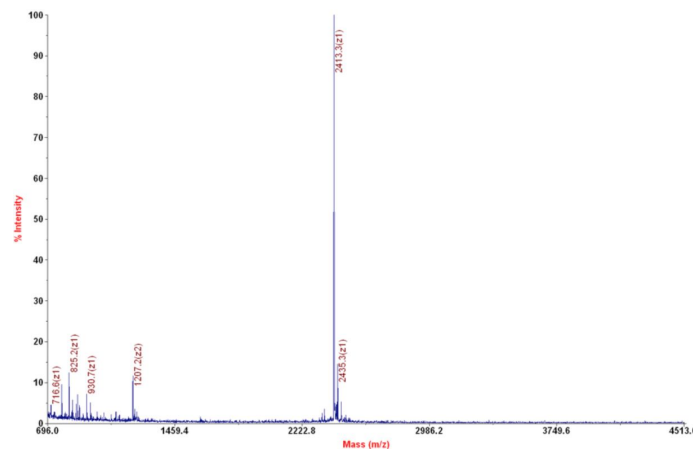
>94% as judged by HPLC and mass spectrometry.

## Storage and Stability:

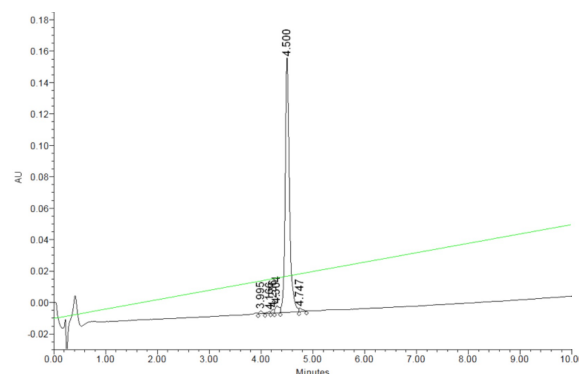
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

Fuchs SM et al (2011). *Curr Biol* 21: 53-58.



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze Histone H2A.ZK4, K8, K12ac Peptide, Biotinylated for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified Histone H2A.ZK4, K8, K12ac Peptide, Biotinylated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.