

# Histone H3 K4,9,14,18Ac Peptide, Biotinylated

Catalog No. 12-0006

Lot No. 12278001

Quantity 50 µg



## Product Description:

A synthetic peptide derived from human histone H3.1, amino acids 1-20, acetylated at lysines 4, 9, 14 and 18, with an added lysine residue containing a biotin moiety at the epsilon-amino group, and a C-terminal amide group (CONH<sub>2</sub>).

## Formulation:

50 µg of lyophilized powder.

## Sequence:

ARTK(Ac)QTARK(Ac)STGGK(Ac)APRK(Ac)QL-K(Biot)-CONH<sub>2</sub>

## Molecular Weight:

2705 Da

## Purity:

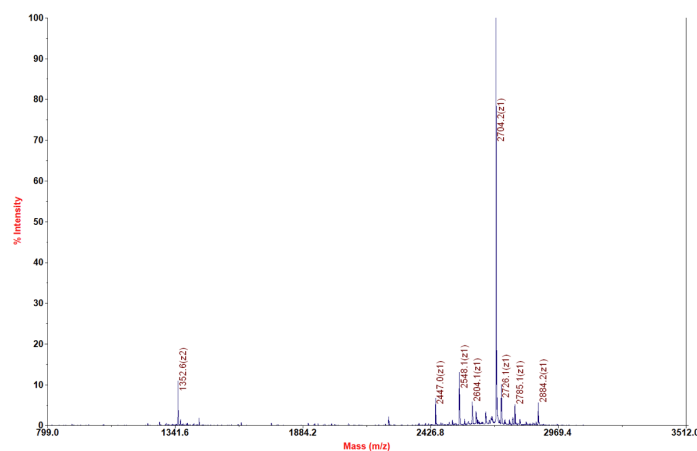
>97% as judged by HPLC and mass spectrometry.

## Storage and Stability:

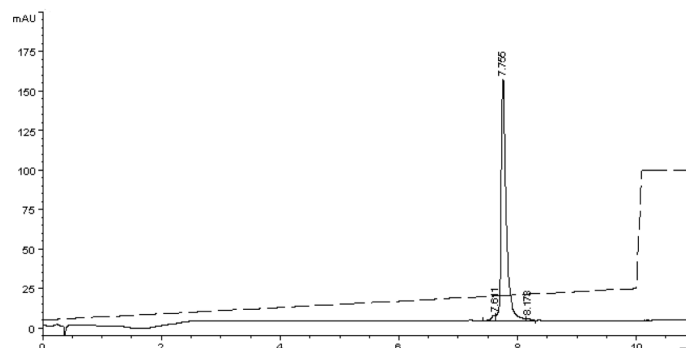
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze histone H3 K4,9,14,18Ac peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H3 K4,9,14,18Ac peptide.