

Histone H4 K5,8,12,16Ac Peptide, Biotinylated

Catalog No. 12-0034

Lot No. 15110002

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H4, amino acids 1-23, acetylated at lysines 5, 8, 12 and 16. The alpha-amino group on the N-terminus is acetylated, with a C-terminal PEG spacer containing a biotin moiety.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-SGRGK(Ac)GGK(Ac)GLGK(Ac)GGAK(Ac)RHRKVLR-Peg (Biot)

Molecular Weight:

2999.5 Da

Purity:

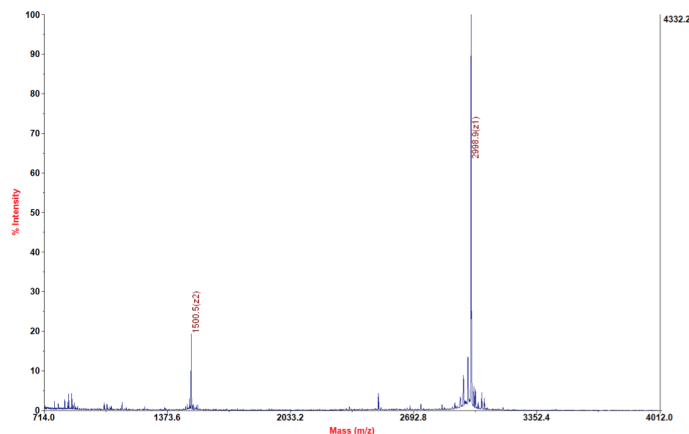
>95.3 % as judged by HPLC and mass spectrometry.

Storage and Stability:

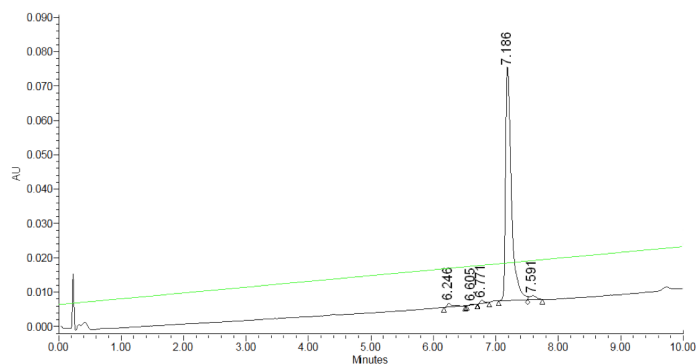
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H4 K5,8,12,16Ac peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H4 K5,8,12,16Ac peptide.