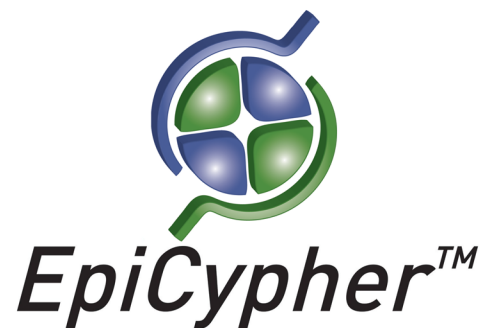


Histone H3 K23Ac Peptide, Biotinylated

Catalog No. 12-0056

Lot No. 14049001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H3.3, amino acids 15-34, acetylated at lysine 23. The alpha-amino group on the N-terminus is acetylated, with a C-terminal PEG spacer containing a biotin moiety at the C-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-APRKQLATK(Ac)AARKSAPSTGG-Peg-Biot

Molecular Weight:

2509 Da

Purity:

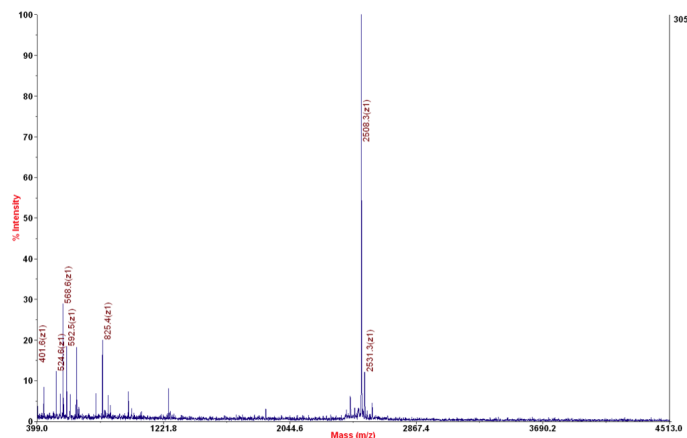
> 97% as judged by HPLC and mass spectrometry.

Storage and Stability:

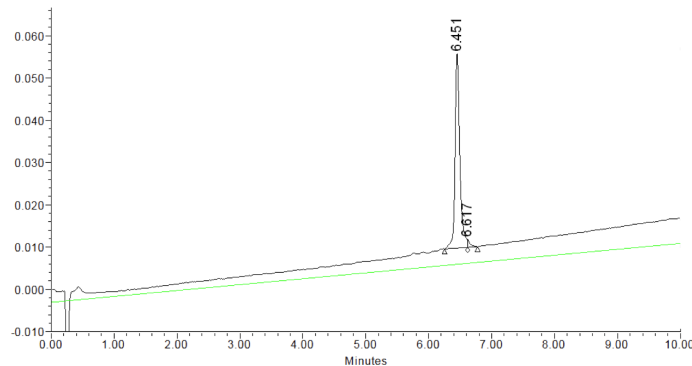
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H3 K23Ac peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H3 K23Ac peptide.