

Histone H4K8, K20ac Peptide, Biotinylated

Catalog No. 12-0135

Lot No. 17251001

Quantity 50 µg



EpiCypher™

Product Description:

A synthetic peptide derived from human histone H4, amino acids 1-23, acetylated at lysines 8 and 20. 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-SGRGKGGK(Ac)GLGKGGAKRHRK(Ac)VLR-Peg(Biot)

Molecular Weight:

2915.5 Da

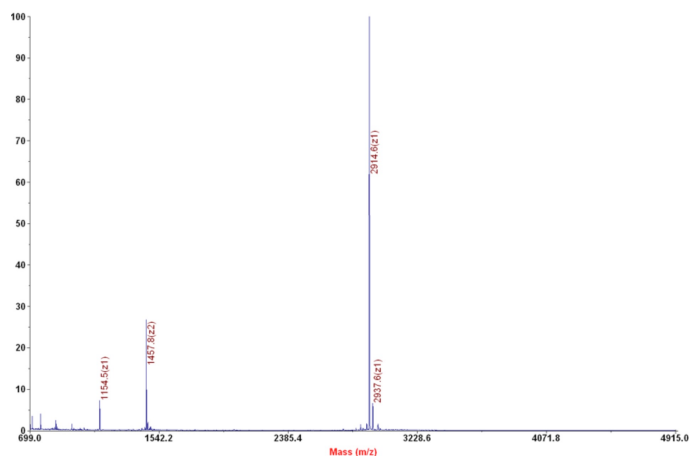
Purity:

>95% as judged by HPLC and mass spectrometry.

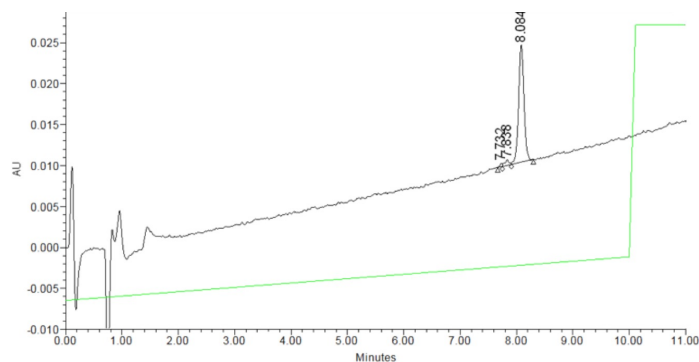
Storage and Stability:

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:



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H4K8, K20ac peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H4K8, K20ac peptide.

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

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com