

# Histone H4K8, K12ac Peptide, Biotinylated

Catalog No. 12-0134

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 12. 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)GLGK(Ac)GGAKRHRKVLR-Peg(Biot)

## Molecular Weight:

2915.5 Da

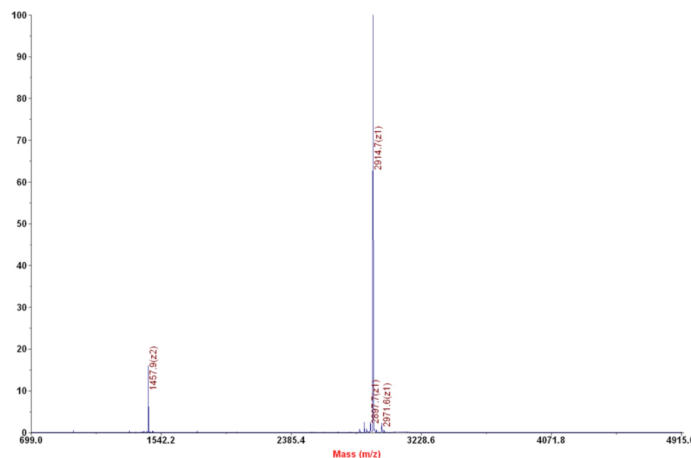
## Purity:

>98% as judged by HPLC and mass spectrometry.

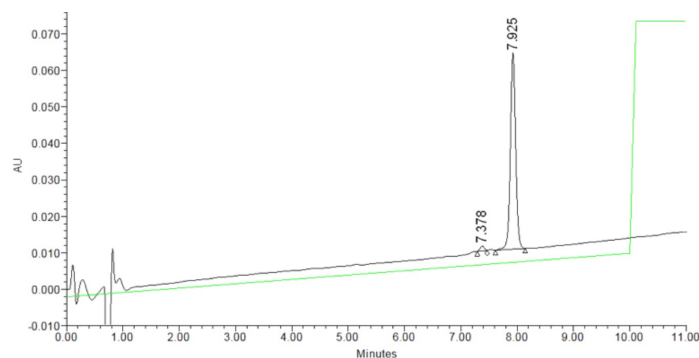
## Storage and Stability:

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, K12ac peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H4K8, K12ac 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](http://www.epicypher.com)