

# Histone H4 N-term Peptide, Biotinylated (No $\alpha$ -Ac)

Catalog No. 12-0057

Lot No. 14224001

Quantity 50  $\mu$ g



## Product Description:

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

## Formulation:

50  $\mu$ g of lyophilized powder.

## Sequence:

SGRGKGGKGLGKGGAKRHRKVLRL-Peg(Biot)

## Molecular Weight:

2789 Da

## Purity:

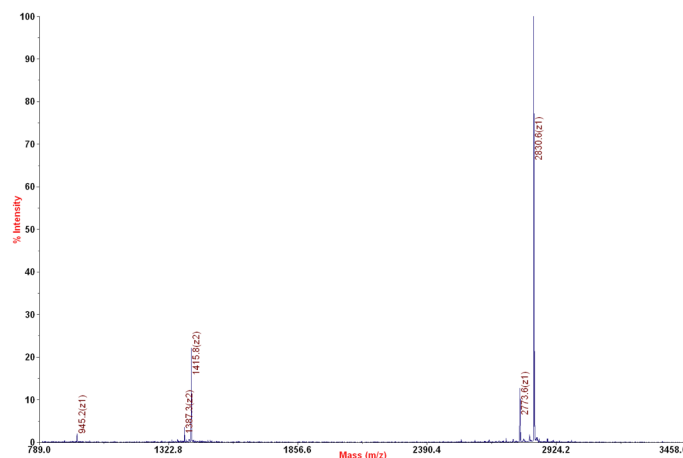
>??% as judged by HPLC and mass spectrometry.

## Storage and Stability:

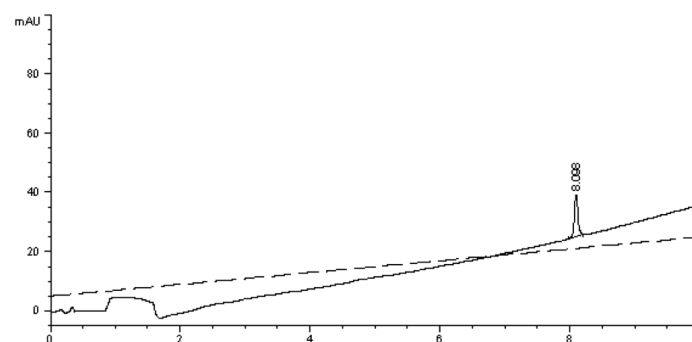
Lyophilized peptides are stable for 2 years at  $-20^{\circ}\text{C}$  from date of shipment. Peptides in solution are stable for 6 months at  $-80^{\circ}\text{C}$ . Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to 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 N-terminal peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H4 N-terminal peptide.