

# Histone H2BK12, K15, K16, K20ac Peptide, Biotinylated

Catalog No. 12-0211

Lot No. 17277001

Quantity 50 µg



EpiCypher™

## Product Description:

A synthetic peptide derived from human histone H2B amino acids 1-24, acetylated at lysines 12, 15, 16, and 20, with a PEG spacer containing a biotin moiety at the C-terminus.

## Formulation:

50 µg of lyophilized powder.

## Sequence:

PEPAKSAPAPKK(Ac)GSK(Ac)K(Ac)AVTK(Ac)AQKK-Peg-Biot

## Molecular Weight:

3072.7 Da

## Purity:

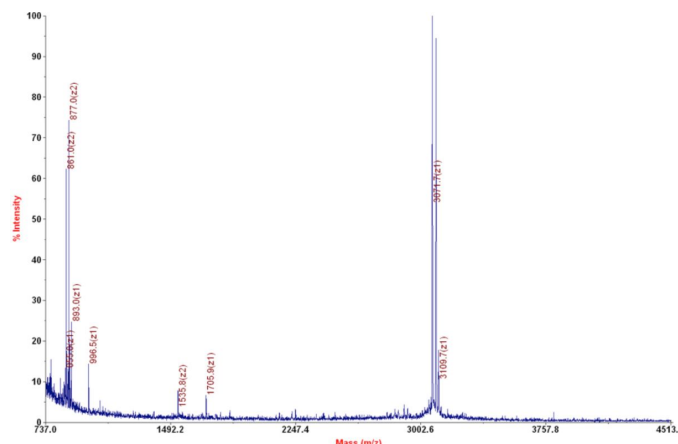
>86% as judged by HPLC and mass spectrometry.

## Storage and Stability:

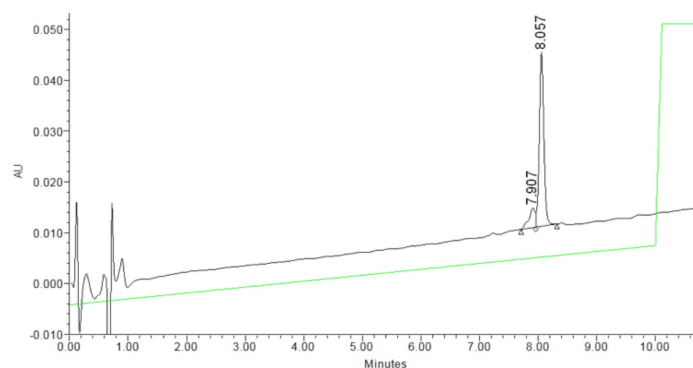
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze Histone H2BK12, K15, K16, K20ac Peptide, Biotinylated for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified Histone H2BK12, K15, K16, K20ac Peptide, Biotinylated.

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