

# Histone H3 aa74-84 Peptide, Biotinylated

Catalog No. 12-0025

Lot No. 12292001

Quantity 50 µg



## Product Description:

A synthetic peptide derived from human histone H3.1, amino acids 74-84. The alpha-amino group on the N-terminus is acetylated, with a PEG spacer containing a biotin moiety and a C-terminal amide group (CONH<sub>2</sub>).

## Formulation:

50 µg of lyophilized powder.

## Sequence:

Ac-IAQDFKTDLRF-Peg(Biot)-CONH<sub>2</sub>

## Molecular Weight:

2051 Da

## Purity:

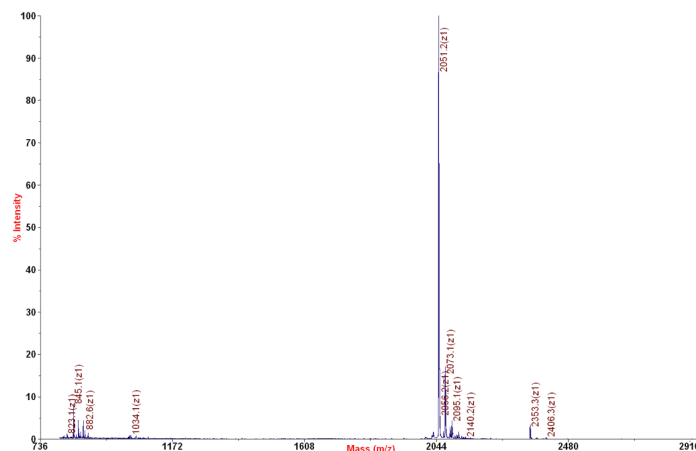
>99% as judged by HPLC and mass spectrometry.

## Storage and Stability:

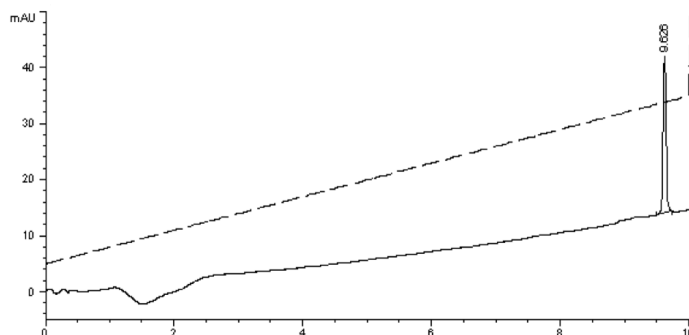
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze histone H3 aa74-84 peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H3 aa74-84 peptide.