

Histone H3 aa27-45 Peptide, Biotinylated

Catalog No. 12-0020

Lot No. 12290001

Quantity 50 µg



Product Description:

A synthetic peptide derived from *S. cerevisiae* histone H3, amino acids 27-45. There are two added glycine residues, an added lysine residue containing a biotin moiety at the epsilon-amino group, and a C-terminal amide group (CONH₂).

Formulation:

50 µg of lyophilized powder.

Sequence:

KSAPSTGGVKKPHRYKPGT-GG-K(Biot)-CONH₂

Molecular Weight:

2464 Da

Purity:

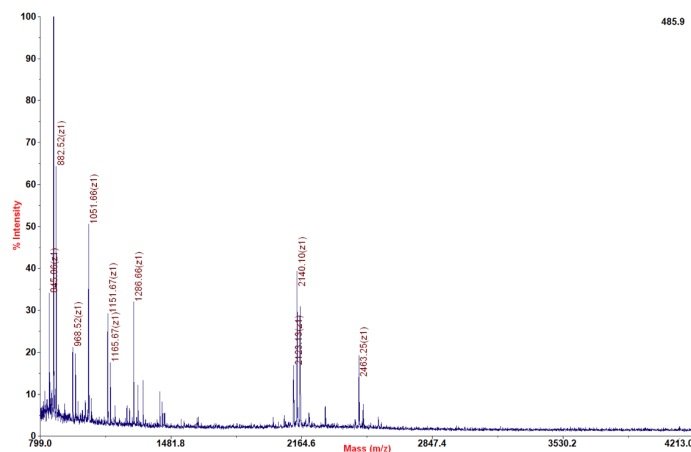
>85% as judged by HPLC and mass spectrometry.

Storage and Stability:

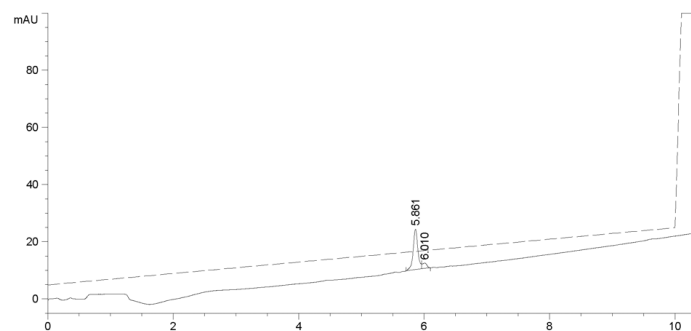
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H3 aa27-45 peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H3 aa27-45 peptide.