

Histone H4 aa11-27 Peptide, Biotinylated

Catalog No. 12-0035

Lot No. 15149002

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H4, amino acids 11-27. 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-GKGGAKRHRKVLRDNIQ-Peg(Biot)

Molecular Weight:

2404 Da

Purity:

>97% 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 H4 aa11-27 peptide for identity and homogeneity.

HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H4 aa11-27 peptide.