

Histone H3K36me3 Peptide - biotinylated

Catalog No: 81057

Lot No: 31017001

Expressed In: Synthetic

Quantity: 100 µg

Source: N/A

Buffer Contents: 100 µg of lyophilized powder.

Background: Active Motif biotinylated histone peptides are suitable for use in a variety of binding, pull-down and enzymatic assays for the study of proteins that interact with histones and that read, write or erase histone modifications.

Protein Details: A synthetic peptide derived from Histone H3.1, amino acids 21-44, with an added lysine residue containing a biotin moiety and a C-terminal amide group (CONH₂). C-term biotinylated Lys residue is separated with two Gly (provide short spacer) from the rest of the peptide sequence. Accurate UV quantification is done based on internal Tyr. The presence of the lysine residue is for quantitation and does not affect peptide functionality.

Peptide Sequence: ATKAARKSAPATGGV-Kme3-KPHRYRPG-GGK(Biotin) - NH₂

Calculated MW Mavg 3033.58

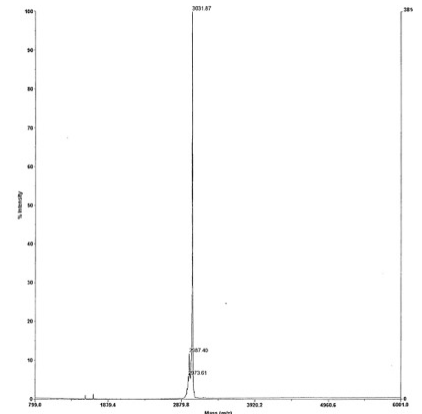
Observed MW Mavg 3031.87

Purity: 95% by HPLC and Mass Spectrometry.

Application Notes: Spin vial briefly in microfuge. Reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot reconstituted peptide to avoid repeated freeze-thawing.

Reference: Fuchs, S.M. *et. al.* Curr Biol. 2011 Jan 11; 21(1): 53–58.

Storage and Guarantee: Store lyophilized peptide at -20°C. Peptides in solution are stable for 6 months at -80°C.



Deconvoluted MALDI-TOF mass spectrum of biotinylated peptide (21-44 H3 histone amino acids).