

Histone H4K20me2 Peptide - biotinylated

Catalog No: 81112

Lot No: 03018001

Expressed In: Synthetic

Buffer Contents: 100 µg of lyophilized powder.

Quantity: 100 µg

Source: N/A

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 H4, amino acids 1-23, 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) and one Tyr (provides accurate UV quantification) from the rest of the peptide sequence. The presence of the tyrosine does not affect peptide functionality.

Peptide Sequence: SGRGKGGKGLGKGGAKRHR-Kme₂-VLRGG-YK(Biotin)-NH₂

Calculated MW Mavg 3019.57

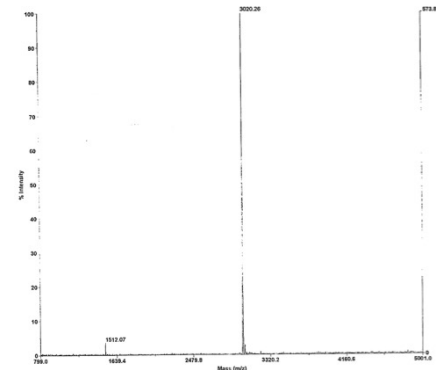
Observed MW Mavg 3020.28

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 (1-23 H4 histone amino acids)