

**Catalog No:** 81038  
**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 1-21, with an added lysine residue containing a biotin moiety and a C-terminal amide group (CONH<sub>2</sub>). 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 lysine residue is for quantitation and does not affect peptide functionality.

Peptide Sequence: ARTKQTARKSTGGKAPRKQLA - GGYK(Biotin) - NH<sub>2</sub>

Calculated MW Mavg 2885.39

Observed MW Mavg 2886.22

Purity: 95% by HPLC and Mass Spectrometry.

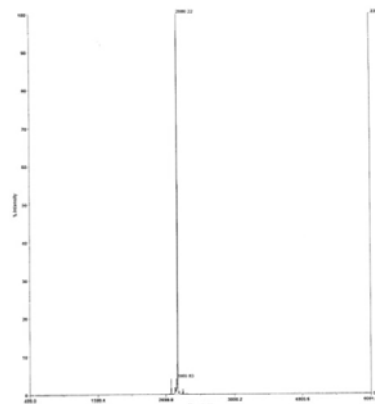
**Application Notes:** Spin vial briefly in microfuge. Reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot reconstituted peptide to avoid repeated freeze-thawing.

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

#### References:

This product was used in the following publications:  
*Mol. Cell.* (2018). 71(1): 25-41. PMID: 29937342.

**Storage and Guarantee:** Store lyophilized peptide at -20°C. Peptides in solution are stable for 6 months at -80°C. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



**Deconvoluted MALDI-TOF mass spectrum of biotinylated peptide (1-21 H3 histone amino acids).**