

# Histone H3K4me3, R8me2a Peptide, Biotinylated

Catalog No. 12-0221

Lot No. 18019001

Quantity 50 µg



EpiCypher™

## Product Description:

A synthetic peptide derived from human histone H3.1, amino acids 1-20, trimethylated at lysine 4 and asymmetrically dimethylated at arginine 8, with an added C-terminal lysine residue containing a biotin moiety.

## Formulation:

50 µg of lyophilized powder.

## Sequence:

ARTK(me3)QTAR(me2a)KSTGGKAPRKQL-K(Biot)-NH<sub>2</sub>

## Molecular Weight:

2608.2 Da

## Purity:

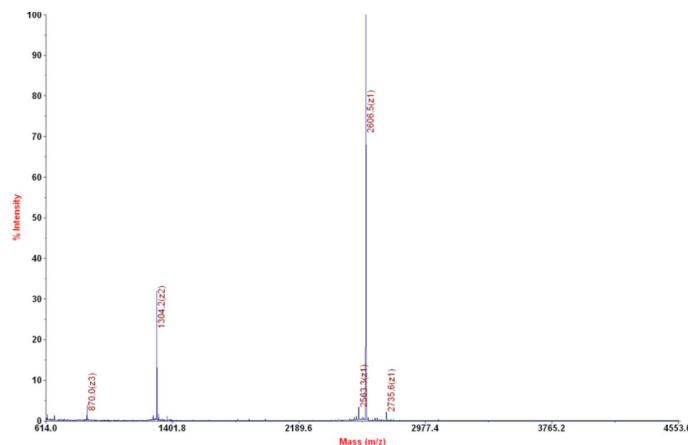
>97% as judged by HPLC and mass spectrometry.

## Storage and Stability:

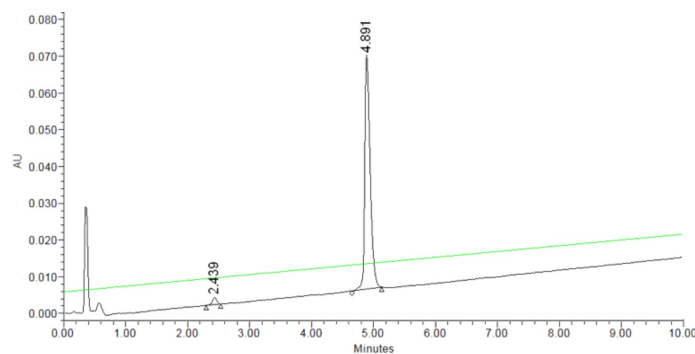
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze histone H3K4me3, R8me2a peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H3K4me3, R8me2a peptide.

This product is for *in vitro* research use only and is not intended for use in humans or animals.