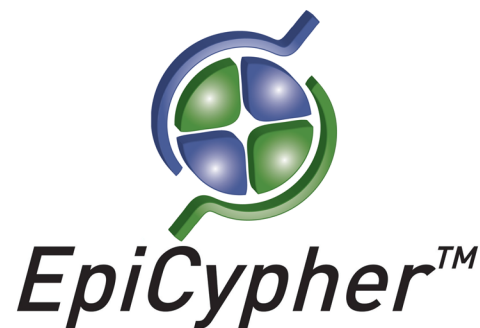


# Histone H3 R2Me2a Peptide, Biotinylated

Catalog No. 12-0060

Lot No. 14224101

Quantity 50 µg



## Product Description:

A synthetic peptide derived from human histone H3, amino acids 1-20, asymmetrically dimethylated at arginine 2. There is an added C-terminal lysine residue containing a biotin moiety at the epsilon amino group, and a C-terminal amide group (CONH<sub>2</sub>)

## Formulation:

50 µg of lyophilized powder.

## Sequence:

AR(Me2a)TKQTARKSTGGKAPRKQL-K(Biot)-CONH<sub>2</sub>

## Molecular Weight:

2565 Da

## Purity:

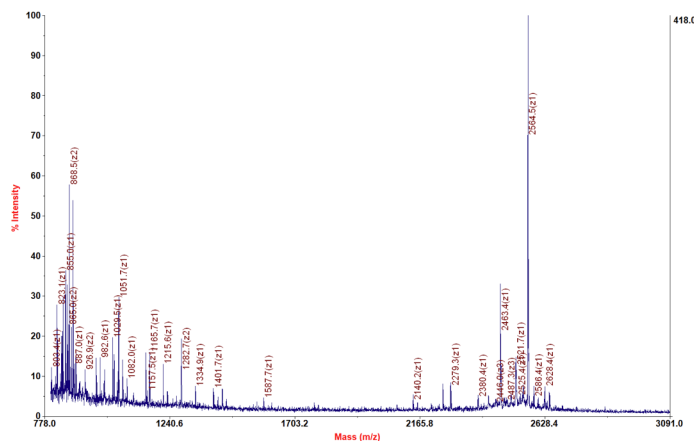
>90% as judged by HPLC and mass spectrometry.

## Storage and Stability:

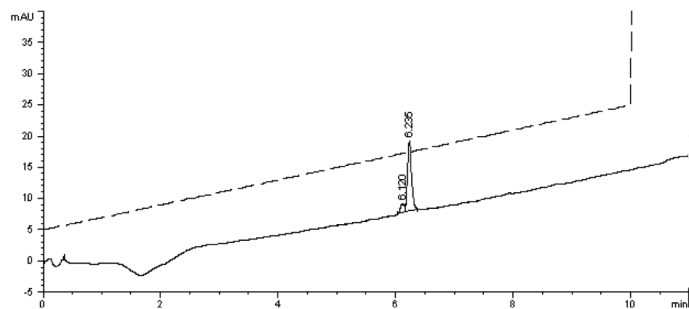
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze histone H3 R2Me2a peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H3 R2Me2a peptide.