

# Histone H4 R3Me2s Peptide, Biotinylated

Catalog No. 12-0059

Lot No. 14224101

Quantity 50 µg



## Product Description:

A synthetic peptide derived from human histone H4, amino acids 1-23, symmetrically dimethylated at arginine 3. The alpha-amino group on the N-terminus is acetylated, and 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:

Ac-SGR(Me2s)GKGGKGLGKGGAKRHRKVLK(Biot)-NH<sub>2</sub>

## Molecular Weight:

2784 Da

## Purity:

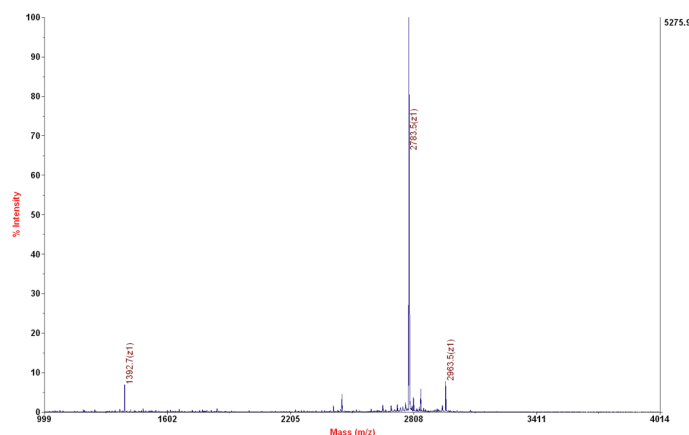
>80% as judged by HPLC and mass spectrometry.

## Storage and Stability:

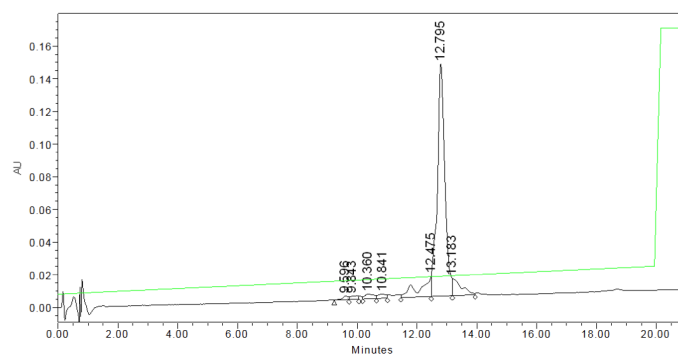
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



**Mass Spec Data:** MALDI-TOF mass spectrometry was used to analyze histone H4 R3Me2s peptide for identity and homogeneity.



**HPLC Data:** Reversed phase analytical HPLC chromatogram of purified histone H4 R3Me2s peptide.