

Histone H4 R3Me2a Peptide, Biotinylated

Catalog No. 12-0058

Lot No. 14224101

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H4, amino acids 1-23, asymmetrically dimethylated on arginine 3. The alpha-amino group on the N-terminus is acetylated, with a C-terminal lysine containing a biotin moiety.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-SGR(Me2a)GKGGKGLGKGGAKRHRKVLR-K(Biot)-NH₂

Molecular Weight:

2784 Da

Purity:

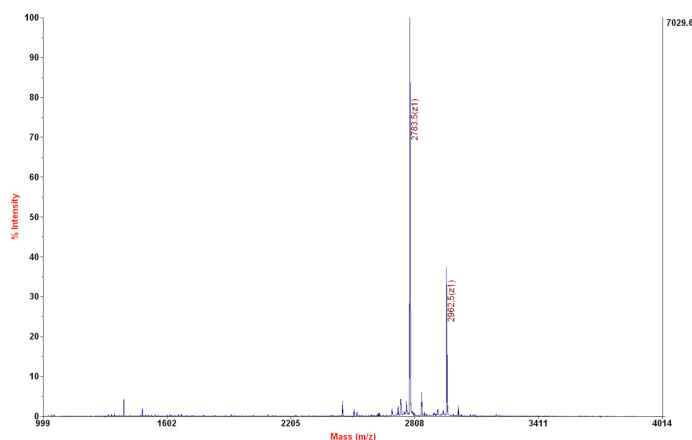
>85% as judged by HPLC and mass spectrometry.

Storage and Stability:

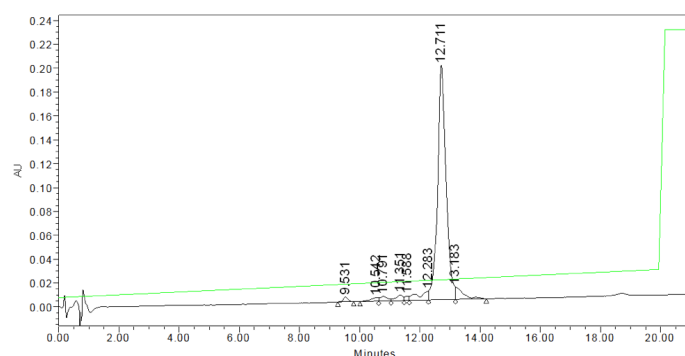
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H4 R3Me2a peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H4 R3Me2a peptide.