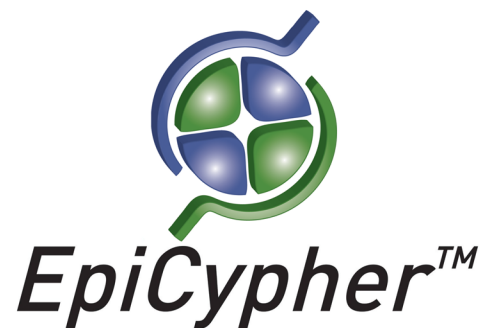


Histone H3 K27Me2 Peptide, Biotinylated

Catalog No. 12-0018

Lot No. 12276001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H3.3, amino acids 15-34, dimethylated at lysine 27. The alpha-amino group on the N-terminus is acetylated, with a PEG spacer containing a biotin moiety at the C-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-APRKQLATKAARK(Me₂)SAPSTGG-Peg(Biot)

Molecular Weight:

2495 Da

Purity:

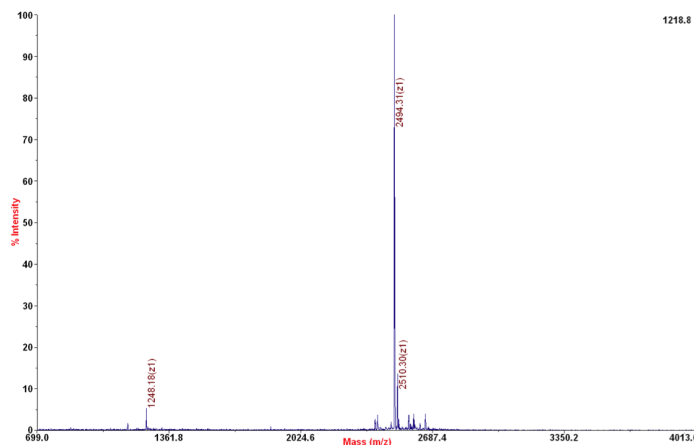
>80% as judged by HPLC and mass spectrometry.

Storage and Stability:

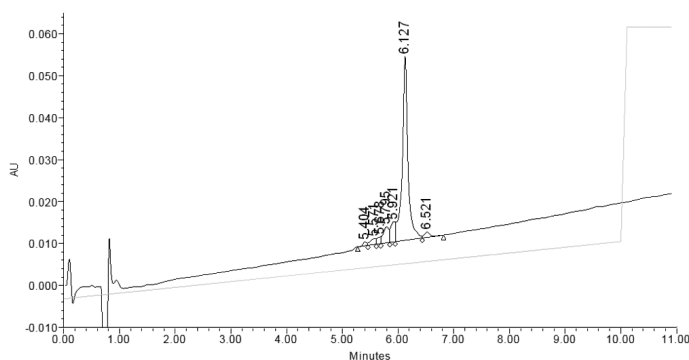
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H3 K27Me2 peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H3 K27Me2 peptide.