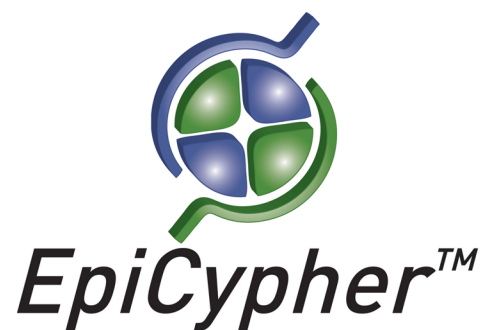


Histone H3 C-terminal Peptide, Biotinylated

Catalog No. 12-0067

Lot No. 15274001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H3.1, amino acids 120-135, with a PEG spacer containing a biotin moiety at the N-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

(Biot)Peg-MPKDIQLARRIGERA-OH

Molecular Weight:

2439 Da

Purity:

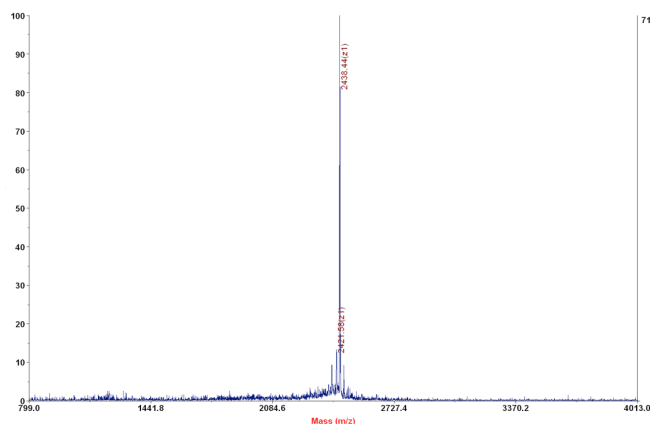
>90% as judged by HPLC and mass spectrometry.

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 C-terminal peptide for identity and homogeneity.