

Histone H3K9cr Peptide, Biotinylated



EpiCypher™

Catalog No. 12-0099
Lot No. 18053001
Pack Size 50 µg

Product Description:

A synthetic peptide derived from human histone H3.1, amino acids 1-20, crotonylated at lysine 9, with an added lysine residue containing a biotin moiety at the epsilon-amino group, and a C-terminal amide group (NH₂).

Formulation:

50 µg of lyophilized powder.

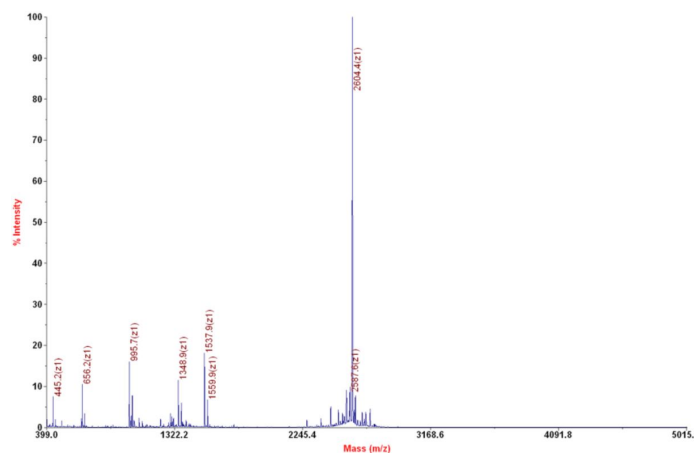
Storage and Stability:

Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6

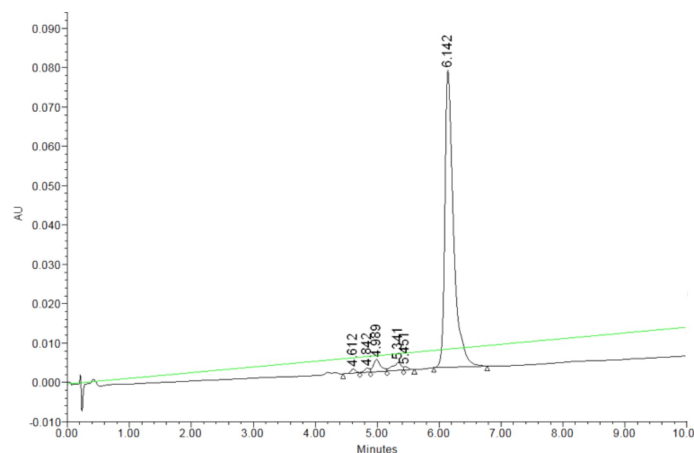
Application Notes:

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H3K9cr peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H3K9cr peptide.