

Histone H2AX N-terminal Peptide, Biotinylated

Catalog No. 12-0062

Lot No. 14283001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone variant H2AX, amino acids 1-17. The N-terminus is acetylated and a PEG spacer containing a biotin moiety on the C-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-SGRGKTGGKARAKAKSR-Peg-Biotin

Molecular Weight:

2187 Da

Purity:

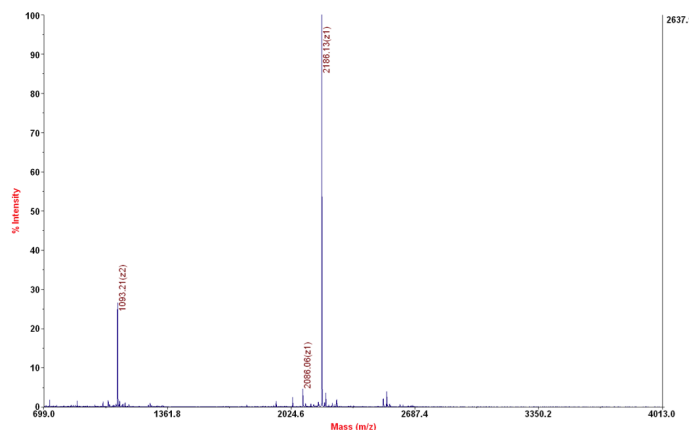
>98% as judged by HPLC and mass spectrometry.

Storage and Stability:

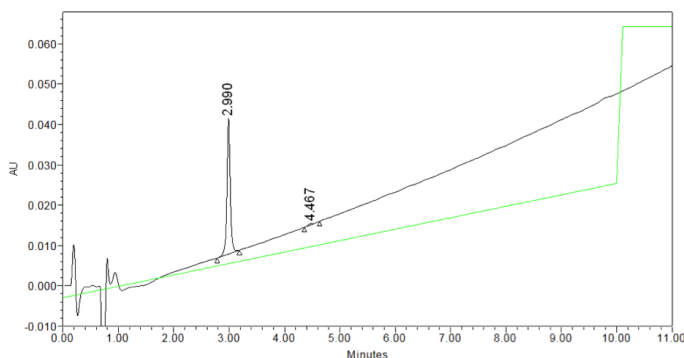
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H2AX N-terminal peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H2AX N-terminal peptide.