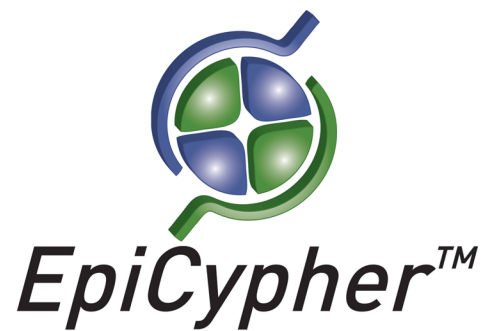


Histone H2AX S139P Peptide, Biotinylated

Catalog No. 12-0040

Lot No. 12289001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone variant H2AX, amino acids 132-142, phosphorylated at serine 139. There is an N-terminal PEG spacer containing a biotin moiety, and a hydroxyl group on the C-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

Peg(Biot)-GKKATQAS(P)QEY-OH

Molecular Weight:

1819 Da

Purity:

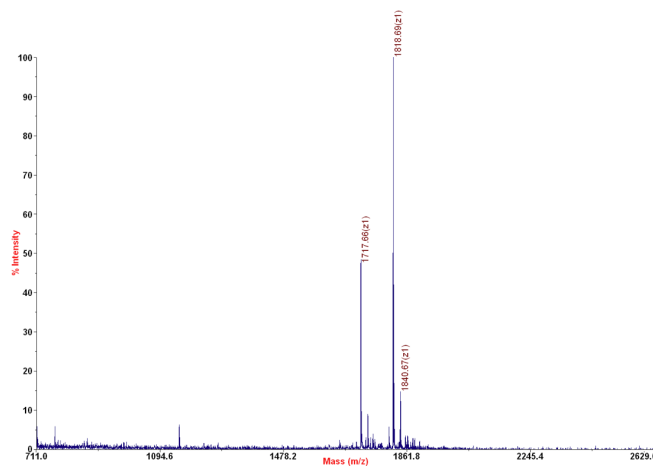
>99% as judged by HPLC and mass spectrometry.

Storage and Stability:

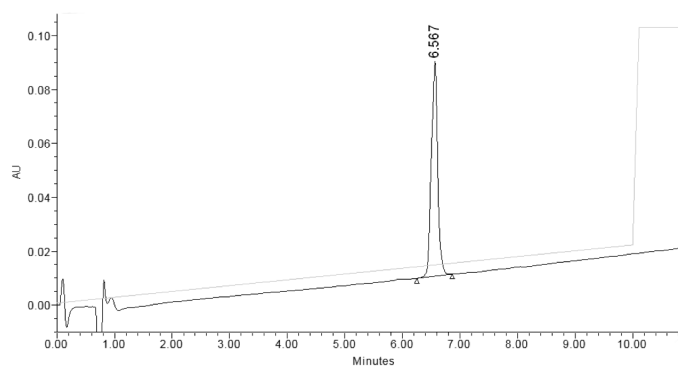
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H2AX S139P peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H2AX S139P peptide.