

# Histone H2AX C-terminal Peptide, Biotinylated

Catalog No. 12-0039

Lot No. 12289001

Quantity 50 µg



## Product Description:

A synthetic peptide derived from human histone variant H2AX, amino acids 132-142. 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)-GKKATQASQEY-OH

## Molecular Weight:

1739 Da

## Purity:

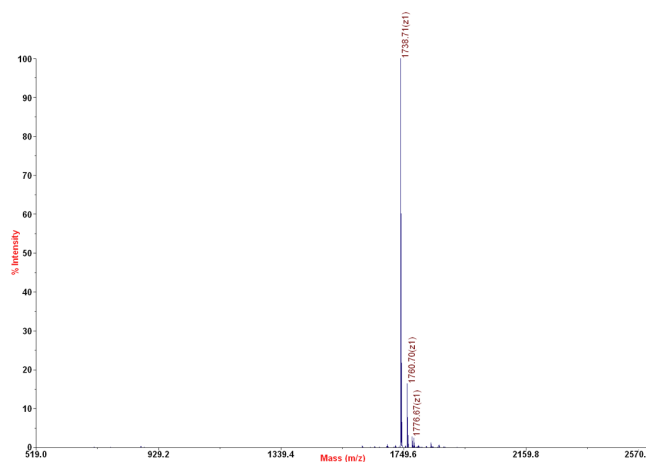
>99% as judged by HPLC and mass spectrometry.

## Storage and Stability:

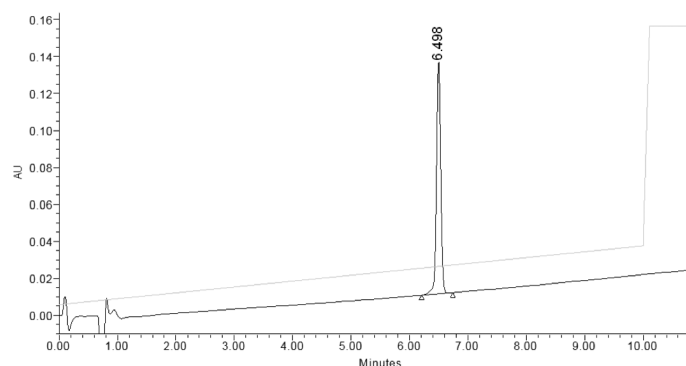
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

## References:

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



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



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