

Histone H3.3T32p, K36me1 Peptide, Biotinylated

Catalog No. 12-0226

Lot No. 18019001

Quantity 50 µg



EpiCypher™

Product Description:

A synthetic peptide derived from human histone H3.3 (accession number P84243), amino acids 27-45, phosphorylated at threonine 32 and monomethylated at lysine 36. There is an added glycine residue, an added lysine residue containing a biotin moiety at the epsilon-amino group.

Formulation:

50 µg of lyophilized powder.

Sequence:

KSAPST(p)GGVK(me1)KPHRYRPGT-G-K(Biot)-NH₂

Molecular Weight:

2528.9 Da

Purity:

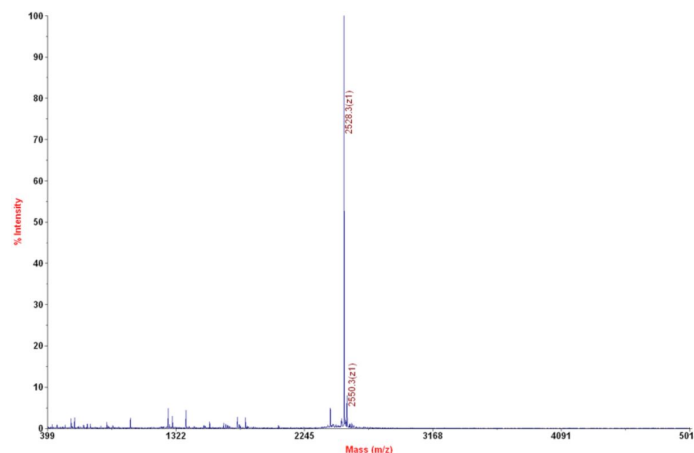
>98% as judged by HPLC and mass spectrometry.

Storage and Stability:

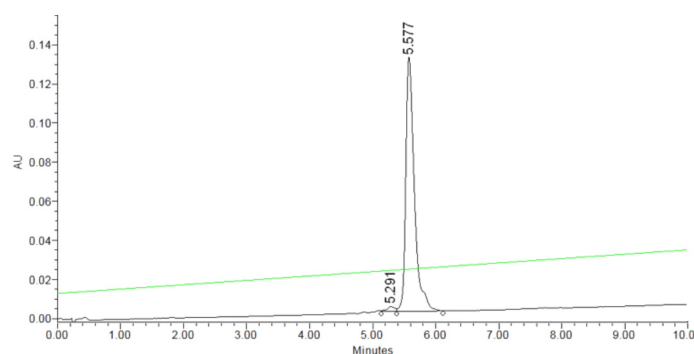
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze histone H3.3T32p, K36me1 peptide for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified histone H3.3T32p, K36me1 peptide.

This product is for *in vitro* research use only and is not intended for use in humans or animals.

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com