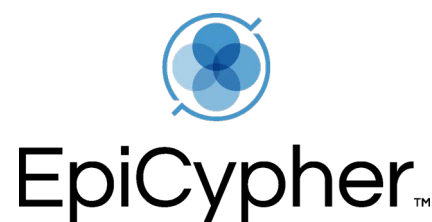


Histone H2AR3me2a, K5ac Peptide, Biotinylated

Catalog No. 12-0192

Lot No. 17277001

Quantity 50 µg



Product Description:

A synthetic peptide derived from human histone H2A amino acids 1-17, asymmetrically dimethylated at arginine 3 and acetylated at lysine 5. The alpha-amino group on the N-terminus is acetylated, with a PEG spacer containing a biotin moiety at the C-terminus.

Formulation:

50 µg of lyophilized powder.

Sequence:

Ac-SGR(Me2a)GK(Ac)QGGKARAKAKTR-Peg-Biot

Molecular Weight:

2297.7 Da

Purity:

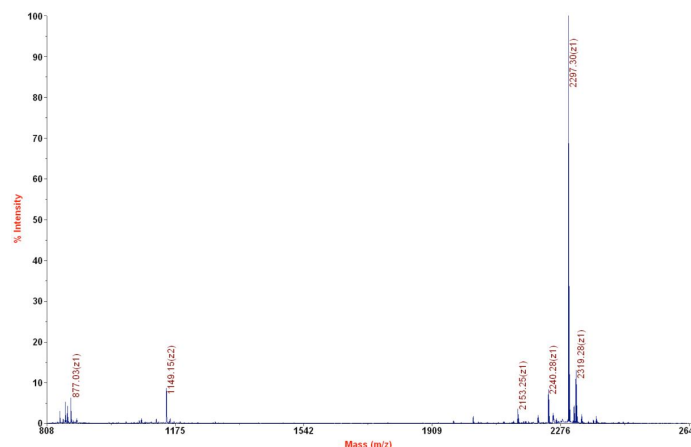
>99% as judged by HPLC and mass spectrometry.

Storage and Stability:

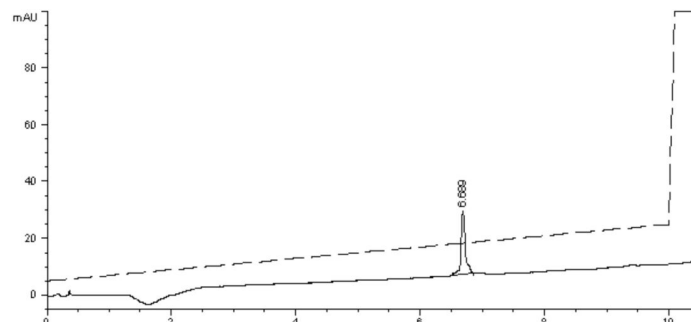
Lyophilized peptides are stable for 2 years at -20°C from date of shipment. Peptides in solution are stable for 6 months at -80°C. Spin vial briefly in microfuge and reconstitute with sterile distilled water and vortex briefly until resuspended. Aliquot peptide to avoid repeated freezing and thawing.

References:

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



Mass Spec Data: MALDI-TOF mass spectrometry was used to analyze Histone H2AR3me2a, K5ac Peptide, Biotinylated for identity and homogeneity.



HPLC Data: Reversed phase analytical HPLC chromatogram of purified Histone H2AR3me2a, K5ac Peptide, Biotinylated.

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