
Histone Acetyl H3K14ac Peptide

(Catalog No. R-1011)

Description: Synthesized acetyl histone H3-K14 peptide derived from residues within 1-100 of human histone H3.

Purity: >90% by HPLC analysis

Concentration: 0.24 mg/ml (200 μ M)

Formulation: PBS buffer

Storage: -20°C , stable for 1 year from the date of shipment. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance.

Application: Enzyme activity
Histone affinity precipitation assay

Research use: Research use only.

Handling Recommendations: For maximum recovery of the products, centrifuge the vial prior to opening the cap.

This product is for research purposes only. Not intended for use in diagnostic procedures.