

Histone Acetyl H3K14ac Peptide, Biotinylated

(Catalog No. R-1012)

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

Purity: >90% by HPLC analysis

Concentration: 0.24 mg/ml (100 µM)

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.

Ordering Information

Products	Size	Cat. No.
Histone Acetyl H3K14ac Peptide, Biotinylated	100 µl	R-1012-100
	200 µl	R-1012-200

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