

Histone Acetyl H3K9ac Peptide

(Catalog No. R-1009)

Description: Synthesized acetyl histone H3-K9 peptide derived from residues within 1-100 of human histone H3.

Purity: >90% by HPLC analysis

Concentration: 0.45 mg/ml (200 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 H3K9ac Peptide	100 µl	R-1009-100
	200 µl	R-1009-200

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