

---

## Histone Acetyl H4K16ac Peptide

(Catalog No. R-1017)

---

**Description:** Synthesized acetyl histone H4-K16 peptide derived from residues within 1-100 of human histone H4.

**Purity:** >90% by HPLC analysis

**Concentration:** 0.4 mg/ml (200  $\mu$ M)

**Formulation:** Distilled water

**Storage:**  $-20^{\circ}\text{C}$ , 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

<b>Products</b>	<b>Size</b>	<b>Cat. No.</b>
Histone Acetyl H4K16ac Peptide	100 $\mu$ l	R-1017-100
	200 $\mu$ l	R-1017-200

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