

---

## Histone Acetyl H4ac Peptide, Biotinylated

(Catalog No. R-1008)

---

|                                  |   |
|----------------------------------|---|
| <b>Description:</b>              | Synthesized acetyl histone H4 (K5, 8, 12, 16) peptide corresponding to within residues 1-18 of human histone H4, biotin conjugated.                       |
| <b>Purity:</b>                   | >90% by HPLC analysis   |
| <b>Concentration:</b>            | 0.25 mg/ml (100 µM)   |
| <b>Formulation:</b>              | Distilled water   |
| <b>Storage:</b>                  | -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. |
| <b>Application:</b>              | Enzyme activity<br>Histone affinity precipitation assay   |
| <b>Research use:</b>             | Research use only.  |
| <b>Handling Recommendations:</b> | 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 H4ac Peptide, Biotinylated | 100 µl      | R-1008-100      |
|   | 200 µl      | R-1008-200      |

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