

---

## Histone H3K79me3 Trimethyl Peptide, Biotinylated

(Cat. No. R-1053)

---

### Description

Synthesized trimethyl histone H3-K79 peptide derived from residues within 1-100 of human histone H3, biotin-conjugated.

### Purity

>90% by HPLC analysis

### Concentration

0.28 mg/ml (100  $\mu$ 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

### 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.*