

---

## Histone H3K9me3 Trimethyl Peptide

(Cat. No. R-1028)

---

<b>Description:</b>	Synthesized trimethyl histone H3-K9 peptide derived from residues within 1-100 of human histone H3.
<b>Purity:</b>	>90% by HPLC analysis
<b>Concentration:</b>	0.54 mg/ml (200 $\mu$ M)
<b>Formulation:</b>	Tris-buffered solution
<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 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 H3K9me3 Trimethyl Peptide	100 $\mu$ l	R-1028-100
	200 $\mu$ l	R-1028-200

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