

---

## Histone H3 Peptide (residues 1-21)

(Cat. No. R-1003)

---

<b>Description:</b>	Synthesized histone H3 peptide corresponding to residues within 1-21 of human histone H3.
<b>Purity:</b>	>90% by HPLC analysis
<b>Concentration:</b>	0.4 mg/ml (200 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>	As substrate for studying enzyme activity and screening inhibitors of histone H3 acetyltransferases (HATs) or methyltransferases (HMTs).
<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 H3 Peptide (residues 1-21)	200 µl	R-1003-200
	400 µl	R-1003-400

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