

## Histone H3 Peptide (residues 21-44), Biotinylated

(Catalog No. R-1006)

<b>Description:</b>	Synthesized histone H3 peptide corresponding to residues within 21-44 of human histone H3, biotin conjugated.
<b>Purity:</b>	>90% by HPLC analysis
<b>Concentration:</b>	0.3 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 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

Products	Size	Cat. No.
Histone H3 Peptide (residues 21-44), Biotinylated	100 µl	R-1006-100
	200 µl	R-1006-200

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