
Histone H3 Peptide (residues 69-89), Biotinylated

(Cat. No. R-1057)

Description:	Synthesized histone H3 peptide corresponding to residues within 69-89 of human histone H3, biotin conjugated.
Purity:	>90% by HPLC analysis
Concentration:	0.28 mg/ml (100 μ M)
Formulation:	Tris-buffer solution
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
Research use:	Research use only.
Handling Recommendations:	For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Ordering Information

Products	Size	Cat. No.
Histone H3 Peptide (residues 69-89), Biotinylated	100 μ l	R-1057-100
	200 μ l	R-1057-200

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