
Biotinylated Phospho Histone Ser 10 Peptide

(Cat. No. R-1047)

Description:	Synthesized phospho histone ser10 peptide derived from residues within 1-100 of human histone H3, biotin conjugated.
Purity:	>90% by HPLC analysis
Concentration:	0.28 mg/ml (100 µM)
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
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.
Biotinylated Phospho Histone Ser 10 Peptide	100 µl	R-1047-100
	200 µl	R-1047-200

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