

---

## Phospho Histone Ser10 Peptide

(Cat. No. R-1046)

---

<b>Description:</b>	Synthesized phospho histone Ser10 peptide derived from residues within 1-100 of human histone H3.
<b>Purity:</b>	>90% by HPLC analysis
<b>Concentration:</b>	0.28 mg/ml (100 µM)
<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>
Phospho Histone Ser10 Peptide	100 µl	R-1046-100
	200 µl	R-1046-200

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