

Histone H3K4me2 Dimethyl Peptide

(Cat. No. R-1020)

Description:	Synthesized dimethyl histone H3-K4 peptide derived from residues within 1-100 of human histone H3.
Purity:	>90% by HPLC analysis
Concentration:	0.54 mg/ml (200 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 H3K4me2 Dimethyl Peptide	100 μ l	R-1020-100
	200 μ l	R-1020-200

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