

<b>PRODUCT NAME</b>			
synthetic peptide: region of H4 containing R3me1			
<b>(Histone H4 methylated at arginine 3)</b>			
<b>Cat. No:</b> sp-993-050	<b>Format:</b> 50 µg	<b>Concentration:</b> 100 pmol/µl	<b>Lot #:</b> 001

**Description:** Synthetic peptide derived from the region of human histone H4 containing the methylated arginine 3.

**MW:** 919.5 kDa

**Format:** In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

**Storage:** Aliquot upon receipt and store at -20°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** April 15, 2010