

<b>PRODUCT NAME</b> synthetic peptide: region of H3 containing the unmodified K4, K9 and K14			
<b>Cat. No:</b> sp-999-050	<b>Format:</b> 50 µg / 50 µl	<b>Concentration:</b> 1 µg/µl	<b>Lot #:</b> 001

**Description:** Synthetic peptide containing the first 16 amino acids of human histone H3 including the unmodified K4, K9 and K14.

**MW:** 1657.9 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