

PRODUCT NAME

synthetic peptide:
region of H4 containing the unmodified K5, K8, K12 and K16

Cat. No: sp-990-050**Format:** 50 µg**Concentration:** 100 pmol/µl**Lot #:** 001

Description: Synthetic peptide containing amino acids 2 to 19 of human histone H4 including the unmodified K5, K8, K12 and K16.

MW: 1777.1 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