

PRODUCT NAME synthetic peptide: region of H3 containing the unmodified K79			
Cat. No: sp-996-050	Format: 50 µg / 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide containing amino acids 71 to 86 of human histone H3 including the unmodified K79.

MW: 1953.2 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