

Synthetic peptide: region of H3 containing the unmodified K4, K9 and K14

Cat. No. C16000999

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 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