

Synthetic peptide: region of H4 containing the unmodified R3 and K5

Cat. No. C16000991

Format: 50 µg

Concentration: C16000991

Lot #: 001

Description: Synthetic peptide containing amino acids 1 to 9 of human histone H4 including the unmodified R3 and K5.

MW: 802.4 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