

Synthetic peptide: region of H4 containing the unmodified R3

(H4R3un)

Cat. No. C16000994

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

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

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.

Availability date: July 31, 2009

Last data sheet update: July 31, 2009

Lot #: 001/ **Synthesis date:** March 26, 2009/ **aliquoting date:** July 31, 2009