

PRODUCT NAME			
synthetic peptide: region of H4 containing K5ac			
(Histone H4 acetylated at lysine 5)			
Cat. No: sp-159-050	Format: 50 µg	Concentration: 100 pmol/µl	Lot #: 001

Description: Synthetic peptide derived from the region of human histone H4 containing the acetylated lysine 5.

MW: 820.4 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