

PRODUCT NAME			
synthetic peptide: region of H3 containing K9/14ac			
(Histone H3 [acetylated lysines 9 and 14])			
Cat. No: sp-005-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide derived from the N-terminal domain of human histone H3 containing the acetylated lysines 9 and 14.

This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K9/14ac (cat# pAb-ACHBHS-044). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K9/14ac: see overview in the Technical Data Sheet of the ChIP-grade Diagenode antibody directed against H3K9/14ac (cat# pAb-ACHBHS-044, also cat# pAb-005-044).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: For long storage, aliquot 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: September 30, 2007

Last data sheet update: September 30, 2007

Lot #: 001/ day of the synthesis: September 11, 2006/ aliquoting day: September 28, 2007