

Synthetic peptide: region of H3 containing K9/14ac

(Histone H3 [acetylated lysines 9 and 14])

Cat. No. C16000005

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 No 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 No pAb-ACHBHS-044, Cat No 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