

PRODUCT NAME			
synthetic peptide: region of H3			
(Histone H3)			
Cat. No: sp-059-050	Format: 50 µg/ 50 µl	Concentration: 1 µg/µl	Lot #: 001

Description: Synthetic peptide derived from the C-terminal domain of human histone H3.

This peptide was used as immunogen to raise the Diagenode antibody directed against histone H3 (cat# pAb-059-050). The same peptide can be used to block this antibody in negative control experiments.

For more information about histone H3: see overview in the Technical Data Sheet of the Diagenode antibody directed against histone H3 (cat# pAb-059-050).

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: May 18, 2007/ aliquoting day: September 28, 2007