

Synthetic peptide: region of H3 containing K18ac

(Histone H3 acetylated at lysine 18)

Cat. No. C16000139

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide corresponding to the region of human histone H3 containing the acetylated lysine 18. This peptide was used as immunogen to raise the Diagenode crude serum directed against H3K18ac (Cat. No. C15310139). The same peptide can be used to block this antibody in negative control experiments. For more information about H3K18ac: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15310139).

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