

Synthetic peptide: region of H4 containing K20me1

(Histone H4 monomethylated at lysine 20)

Cat. No. C16000034

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide derived from the region of human histone H4 containing the monomethylated lysine 20. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H4K20me1 (Cat. No. C15310034). The same peptide can be used to block this antibody in negative control experiments.

For more information about H4K20me1: see the Technical Data Sheet of the corresponding Diagenode antibody.

MW: 1638 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 14, 2010