

Synthetic peptide: region of H3 containing K27me2

(Histone H3 dimethylated at lysine 27)

Cat. No. C16000046

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide derived from the region of human histone H3 containing the dimethylated lysine 27. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K27me2 (Cat. No. C15410046). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K27me2: see the Technical Data Sheet of the corresponding Diagenode antibody.

MW: 1417.7 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