

## Synthetic peptide: region of H3 containing K27me1

(Histone H3 dimethylated at lysine 27)

**Cat. No. C16000045**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide derived from the region of human histone H3 containing the monomethylated lysine 27. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K27me1 (Cat. No. C15410045). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K27me1: see the Technical Data Sheet of the corresponding Diagenode antibody.

**MW:** 1403.7 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