

## Synthetic peptide: region of H3 containing K4me3

(Histone H3 trimethylated at lysine 4)

**Cat. No. C16000003**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide derived from the region of human histone H3 containing the trimethylated lysine 4. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K4me3 (Cat. No. C15410003, C15410203, C15310003). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K4me3: see the Technical Data Sheet of the corresponding Diagenode antibody.

**MW:** 1076.2 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 16, 2010