

## Synthetic peptide: region of H4 containing K20me2

(Histone H4 dimethylated at lysine 20)

Cat. No. **C16000138**

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

**Description:** Synthetic peptide derived from the region of human histone H4 containing the dimethylated lysine 20

**MW:** 1652 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 15, 2010