

## **Synthetic peptide: region of H3 containing the unmodified K23, K27 and S28**

**Cat. No.** C16000998

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide containing amino acids 20 to 35 of human histone H3 including the unmodified K23, K27 and S28.

**MW:** 1498.8 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