

## **Synthetic peptide: region of H4 containing the unmodified R3me1**

(Histone H4 methylated at arginine 3)

**Cat. No.** C16000993

**Format:** 50 µg

**Concentration:** 100 pmol/µl

**Lot #:** 001

**Description:** Synthetic peptide derived from the region of human histone H4 containing the methylated arginine 3.

**MW:** 919.5 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