

## Synthetic peptide: region of H4 containing K12/16/20ac

(Histone H4 acetylated at lysine 12, 16 and 20)

**Cat. No. C16000164**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide derived from the region of human histone H4 containing the acetylated lysines 12, 16 and 20.

**MW:** 1266.4 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