

**PRODUCT NAME**

synthetic peptide:  
region of H4 containing K12ac

(Histone H4 acetylated at lysine 12)

**Cat. No:** sp-162-050

**Format:** 50 µg

**Concentration:** 100 pmol/µl

**Lot #:** 001

**Description:** Synthetic peptide derived from the region of human histone H4 containing the acetylated lysine 12.

**MW:** 762.5 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