

## Synthetic peptide: region of H4 containing R3me2(asym)

(Histone H4 asymmetrically dimethylated at arginine 3)

**Cat. No.** C16000150

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

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

**MW:** 935.2 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