

|   |                              |                               |                   |
|---|------------------------------|-------------------------------|-------------------|
| <b>PRODUCT NAME</b>                                 |                              |                               |                   |
| synthetic peptide:<br>region of H3 containing K14ac |                              |                               |                   |
| [Histone H3 acetylated at lysine 14]                |                              |                               |                   |
| <b>Cat. No:</b> sp-131-050                          | <b>Format:</b> 50 µg / 50 µl | <b>Concentration:</b> 1 µg/µl | <b>Lot #:</b> 001 |

**Description:** Synthetic peptide derived from the region of human histone H3 containing the acetylated lysine 14.

**MW:** 918 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