

| | | | |
|---|------------------------------|-------------------------------|-------------------|
| PRODUCT NAME synthetic peptide: region of H4 containing K20me2 | | | |
| (Histone H4 dimethylated at lysine 20) | | | |
| Cat. No: sp-138-050 | Format: 50 µg / 50 µl | Concentration: 1 µg/µl | Lot #: 001 |

Description: Synthetic peptide derived from the region of human histone H4 containing the dimethylated lysine 20

MW: 1652 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