

| | | | |
|--|------------------------------|-------------------------------|-------------------|
| PRODUCT NAME | | | |
| synthetic peptide: region of H3 containing K36me3 | | | |
| (Histone H3 trimethylated at lysine 36) | | | |
| Cat. No: sp-058-050 | Format: 50 µg / 50 µl | Concentration: 1 µg/µl | Lot #: 001 |

Description: Synthetic peptide derived from the region of human histone H3 containing the trimethylated lysine 36. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K36me3 (cat# pAb-058-050, CS-058-100). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K36me3: see the Technical Data Sheet of the corresponding Diagenode antibody.

MW: 1486.7 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 16, 2010