
Histone H3K36me1 Monomethyl Peptide, Biotinylated

(Cat. No. R-1037)

| | |
|----------------------------------|---|
| Description: | Synthesized monomethyl histone H3-K36 peptide derived from residues within 1-100 of human histone H3, biotin conjugated. |
| Purity: | >90% by HPLC analysis |
| Concentration: | 0.3 mg/ml (100 µM) |
| Formulation: | PBS buffer |
| Storage: | -20°C, stable for 1 year from the date of shipment. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. |
| Application: | Enzyme activity Histone affinity precipitation assay |
| Research use: | Research use only. |
| Handling Recommendations: | For maximum recovery of the products, centrifuge the vial prior to opening the cap. |

Ordering Information

| Products | Size | Cat. No. |
|---|--------|------------|
| Histone H3K36me1 Monomethyl Peptide, Biotinylated | 100 µl | R-1037-100 |
| | 200 µl | R-1037-200 |

This product is for research purposes only. Not intended for use in diagnostic procedures.