
Histone H4K20me2 Dimethyl Peptide, Biotinylated

(Cat. No. R-1055)

Description

Synthesized dimethyl histone H4-K20 peptide derived from residues within 1-100 of human histone H4, biotin-conjugated.

Purity

>90% by HPLC analysis

Concentration

0.28 mg/ml (100 μ M)

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

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

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