

## **Product Data Sheet**

Product Name: [Lys(Me<sub>3</sub>)<sup>9</sup>] - Histone H3 (1 - 21) - GGK(Biotin), H3K9(Me<sub>3</sub>), biotin-labeled

Catalog Number: AS-64360-025 (0.25 mg) Lot Number: See label on vial

AS-64360-1 (1 mg)

Sequence: H-Ala-Arg-Thr-Lys-Gln-Thr-Ala-Arg-Lys(Me3)-Ser-Thr-Gly-Gly-Lys-Ala-

Pro-Arg-Lys-Gln-Leu-Ala-Gly-Gly-Lys(Biotin)-OH (3-letter codse)
ARTKQTAR-K(Me3)-STGGKAPRKQLA-GGK(Biotin) (1-letter code)

Molecular Weight: 2765.2 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at – 20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: This peptide is Histone H3 amino acid residues 1 to 21 tri-methylated at Lys-9 with a C-terminal GG linker followed by a biotinylated lysine. Provided at >95% peptide purity, this peptide was dissolved in distilled water at 1 mg/ml and re-lyophilized to powder form.

## **Related Products:**

Click on the following links for a complete Histone listings:

Histone H1, H2A & H2B  $GO^{\otimes}$  Peptides Histone H3 (1-21)  $GO^{\otimes}$  Peptides Histone H3 (21-44)  $GO^{\otimes}$  Peptides Histone H4  $GO^{\otimes}$  Peptides

For Research Use Only