



Product Data Sheet

Product Name: [Lys(Me₂)⁹]-Histone H3 (1-21)-GGK(Biotin), H3K9(Me₂), biotin-labeled
Catalog Number: AS-64359-025 (0.25 mg) Lot Number: See label on vial
AS-64359-1 (1 mg)
Sequence: H-Ala-Arg-Thr-Lys-Gln-Thr-Ala-Arg-Lys(Me₂)-Ser-Thr-Gly-Gly-Lys-Ala-
Pro-Arg-Lys-Gln-Leu-Ala-Gly-Gly-Lys(Biotin)-OH (3-letter code)
ARTKQTAR-K(Me₂)-STGGKAPRKQLA-GGK(Biotin) (1-letter code)
Molecular Weight: 2751.2

% 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 di-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[®] Peptides](#)
[Histone H3 \(1-21\) GO[®] Peptides](#)
[Histone H3 \(21-44\) GO[®] Peptides](#)
[Histone H4 GO[®] Peptides](#)

For Research Use Only