



## Product Data Sheet

**Product Name:** [Lys(Me<sub>2</sub>)<sup>4</sup>]-Histone H3 (1-21)-GGK(Biotin), H3K4(Me<sub>2</sub>), biotin-labeled  
**Catalog Number:** AS-64356-025 (0.25 mg) Lot Number: See label on vial  
AS-64356-1 (1 mg)  
**Sequence:** H-Ala-Arg-Thr-Lys(Me<sub>2</sub>)-Gln-Thr-Ala-Arg-Lys-Ser-Thr-Gly-Gly-Lys-Ala-  
Pro-Arg-Lys-Gln-Leu-Ala-Gly-Gly-Lys(Biotin)-OH (3-letter sequence)  
ART-K(Me<sub>2</sub>)-QTARKSTGGKAPRKQLA-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-4 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<sup>®</sup> Peptides](#)  
[Histone H3 \(1-21\) GO<sup>®</sup> Peptides](#)  
[Histone H3 \(21-44\) GO<sup>®</sup> Peptides](#)  
[Histone H4 GO<sup>®</sup> Peptides](#)

***For Research Use Only***