



## Product Data Sheet

---

**Product Name:** [Lys(Me<sub>3</sub>)<sup>27</sup>]-Histone H3 (21-44)-GK(Biotin), H3K27(Me<sub>3</sub>), biotin-labeled

**Catalog Number:** 64367-025 (0.25 mg)      Lot Number: See label on vial  
64367-1 (1 mg)

**Sequence:** H-Ala-Thr-Lys-Ala-Ala-Arg-Lys(Me<sub>3</sub>)-Ser-Ala-Pro-Ala-Thr-Gly-Gly-Val-Lys-Lys-Pro-His-Arg-Tyr-Arg-Pro-Gly-Gly-Lys(Biotin)-OH (3-letter sequence)  
ATKAAR-K(Me<sub>3</sub>)-SAPATGGVKKPHRYRPG-GK(Biotin) (1-letter code)

**Molecular Weight:** 2959.5

**% Peak Area by HPLC:** ≥ 95

**Appearance:**

**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.

**Related Products:**

Click on the following links for 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***