



## Product Data Sheet

**Product Name:** Histone H3 (1-21), Biotinylated  
**Catalog Number:** AS-61702 (1 mg) **Lot Number:** See label on vial  
**Sequence:** Ala-Arg-Thr-Lys-Gln-Thr-Ala-Arg-Lys-Ser-Thr-Gly-Gly-Lys-Ala-Pro-Arg-Lys-Gln-Leu-Ala-Gly-Gly-Lys(Biotin)-NH<sub>2</sub> (3-letter code)  
ARTKQTARKSTGGKAPRKQLA-GG-K(BIOTIN)-NH<sub>2</sub> (1-letter code)

**Molecular Weight:** 2722.2

**% 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 sequence corresponds to the first 21 amino acids of the NH<sub>2</sub> terminal of histone H3 followed by a GG linker and a biotinylated lysine. This peptide was used to investigate the characteristics and mechanisms of ethanol-induced histone H3 acetylation in rat hepatocytes. Immunocytochemical and immunoblot analyses revealed that ethanol treatment significantly increased H3 acetylation at Lys9 with negligible effects at Lys14, 18, and 23. Ref: Park, P. et al. *Am J Physiol Gastrointest Liver Physiol* **289**, G1124 (2005).

**Additional Information:** Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Biotinylated histone peptides were purchased from AnaSpec. Five micrograms of purified proteins were incubated with 0.2 µg of biotinylated histone peptide in 100 µl of the binding buffer (50 mM Tris-HCl [pH7.5], 150mM NaCl, 0.05% NP-40, 0.3 mg/ml BSA) overnight at 4°C-[Iwase, S. et al. J Cell 128, 1077 \(2007\).](#)

**Published Citations:**

[Iwase, S. et al. J Cell 128, 1077 \(2007\).](#)

**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**