



## Product Data Sheet

---

**Product Name:** Histone H3 (1-21), N-Terminal  
**Catalog Number:** AS-61701 (1 mg) **Lot Number:** See label on vial  
**Sequence:** H-Ala-Arg-Thr-Lys-Gln-Thr-Ala-Arg-Lys-Ser-Thr-Gly-Gly-Lys-Ala-Pro-Arg-Lys-Gln-Leu-Ala-OH (3-letter sequence)  
ARTKQTARKSTGGKAPRKQLA (1-letter code)  
**Molecular Weight:** 2254.6

**% 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.

### 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***