



Product Data Sheet

Product Name: Histone H3 (1-25), amide

Catalog Number: AS-61703 (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-Thr-Lys-Ala-Ala-NH₂ (3-letter code)
ARTKQTARKSTGGKAPRKQLATKAA-NH₂ (1-letter code)

Molecular Weight: 2625.1

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Histone H3 (1-25), amide peptide is freely soluble in water.

Storage: Histone H3 (1-25), amide peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: This sequence corresponds to the N-terminus of histone H3, spanning amino acids 1 to 25, amidated. Histones are small proteins (11–22 kDa) that mediate the folding of DNA into chromatin. Ref: Kobza, K. et al. FEBS J. **272**, 4249 (2005).

Related Products:

Name	Cat #	Size
Histone H3 (1-21), N-Terminal ARTKQTARKSTGGKAPRKQLA	AS-61701	1 mg

For Research Use Only