



## Product Data Sheet

---

<b>Product Name:</b>	Histone H3 (1-20), N-Terminal	
<b>Catalog Number:</b>	AS-62753 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Ala-Arg-Thr-Lys-Gln-Thr-Ala-Arg-Lys-Ser-Thr-Gly-Gly-Lys-Ala-Pro-Arg-Lys-Gln-Leu-OH (3-letter code) ARTKQTARKSTGGKAPRKQL-NH2 (1-letter code)	
<b>Molecular Weight:</b>	2183.6	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Histone H3 (1-20) peptide is freely soluble in water.

**Storage:** Histone H3 (1-20) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This is amino acids 1 to 20 fragment of the histone H3. Comparison the acetylation efficiency of different substrates showed that this peptide corresponding to the N-terminal of H3 histone has nearly identical acetylation efficiency as the H4 peptides. Acetylation of histones is generally associated with active transcription, constitutes a post-translational mark recognized by specific chromatin factors, and has been shown in vitro to prevent salt-induced folding of nucleosome arrays. Multisubunit histone acetyltransferase (HAT) complexes recognize and perform efficient acetylation on nucleosome substrates. Ref: Berndsen, C. et al. *Biochem.* **46**, 2091 (2007).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
[Lys(Ac)9]-Histone H3 (1-20), H3K9(Ac) ARTKQTAR-K(Ac)-STGGKAPRKQL	AS-63680	1 mg
Histone H3 (1-21), N-Terminal ARTKQTARKSTGGKAPRKQLA	AS-61701	1 mg
Histone H3 (1-21), Biotinylated ARTKQTARKSTGGKAPRKQLA-GG-K(BIOTIN)-NH2	AS-61702	1 mg
Histone H3 (1-21), FAM labeled ARTKQTARKSTGGKAPRKQLAGG-K(FAM)-NH2	AS-63824	1 mg

Histone H3 (1-25), amide  
ARTKQTARKSTGGKAPRKQLATKAA-NH<sub>2</sub>

AS-61703 1 mg

*For Research Use Only*