



Product Data Sheet

Product Name: Histone H4 (1-21), p300/CBP Substrate

Catalog Number: AS-62499 (1 mg) **Lot Number:** See label on vial

Sequence: H-Ser-Gly-Arg-Gly-Lys-Gly-Gly-Lys-Gly-Leu-Gly-Lys-Gly-Gly-Ala-Lys-Arg-His-Arg-Lys-Val-OH (3-letter code)
SGRGKGGKGLGKGGAKRHRKV (1-letter code)

Molecular Weight: 2091.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Peptide YY is freely soluble in water.

Storage: Peptide YY 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 is amino acids 1 to 21 fragment of the histone 4. Histone acetyltransferase (HAT) p300/CBP catalyses the acetylation of this N-terminal tail of free histone H4 at Lys5, 8, 12, 16. Lys8 is thought to be the preferred acetylation site in H4. Determinants of interaction of p300 with histone 4 appear to be fully present on H4-21, the N-terminal region of the protein. Ref: Sun, Y. et al. *Mol. Cell. Biol.* Oct 2007, 10.1128/MCB.01382-07.

Related Products:

Name	Cat #	Size
Histone H4 (1-21)-Lys(Biotin)	AS-62555	1 mg
Ac-SGRGKGGKGLGKGGAKRHRKVGG-K(Biotin)		

For Research Use Only