

Product Information

Description:	Human Histone 4 (HIST2H4A), full length encompassing amino acids 2-103(end). The construct contains an N-terminal His-tag (6xHis). This recombinant protein was affinity purified.
Background:	Histones are nuclear proteins forming the nucleosome structure of the chromosomes in eukaryotes, which consists of DNA (Deoxyribonucleic acid) wrapped around an octamer composed of two each of the four core histones H2A, H2B, H3, and H4. The chromatin fiber is further compacted through interaction of histone H1. Histones are modified post-transcriptionally by methylation and acetylation to regulate gene transcription. Purified histones are used as substrates to study the role of histone-modifying enzymes including methyltransferases, acetyltransferases, deacetylases, and demethylases, involved in epigenetic programming.
Species:	Human
Construct:	Histone H4 (His-2-103(end))
Concentration:	0.64 mg/ml
Expression System:	<i>E. coli</i>
Purity:	75%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, and 20% glycerol
MW:	12 kDa
Genbank Accession:	NM_003548
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful as substrate for histone methyltransferase and acetyltransferase assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

