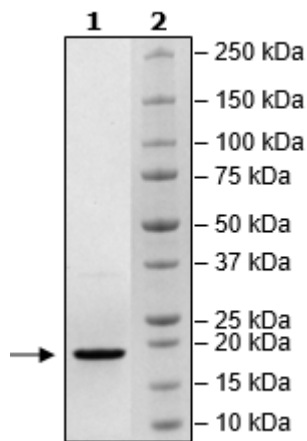


Product Information

Construct:	Histone H3 (C97A, C111A) (MAC-His-2-136(end))-(Biotin)
Mutations:	C97A, C111A
Label:	Histone H3 was modified to contain an N-terminal Met-Ala-Cys, and two additional mutations (C97A, C111A) removed the protein cysteine residues. This ensures site-specific biotinylation at the N-terminal cysteine. Biotin was added chemically to this free cysteine.
Concentration:	0.74 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_003532
MW:	16 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

