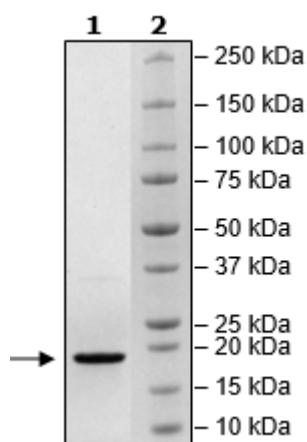


## Product Information

<b>Construct:</b>	Histone H3 (C97A, C111A) (MAC-His-2-136(end))-(Biotin)
<b>Mutations:</b>	C97A, C111A
<b>Label:</b>	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.
<b>Concentration:</b>	0.74 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_003532
<b>MW:</b>	16 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

