

Product Information

Construct:	Histone H2a (His-2-130(end))
Concentration:	5.31 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 8.0, 150 mM NaCl, 10% glycerol, 1 mM PMSF and 1 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_033445
MW:	15 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

