

Product Information

Construct:	Histone H1 (His-2-194(end))
Concentration:	2.12 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, and 20% glycerol.
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_005318
MW:	22 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

