

Product Information

Construct:	Human core histones (two copies each of H2A, H2B, H3, and H4), purified from cultured HeLa cells. Histones are separated from DNA under high salt conditions.
Concentration:	1.30 mg/ml
Species:	Human
Formulated In:	50 mM sodium phosphate buffer, pH 6.8, 2.5 M NaCl
Expression System:	HeLa
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Global MW:	110 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

