

H2A-H2B dimer

CATALOG NO.: HMT-12-316

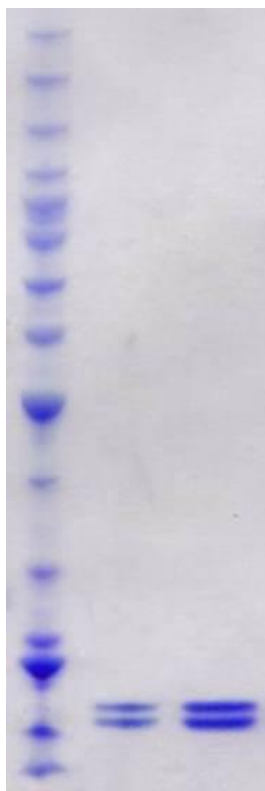
LOT NO.:

DESCRIPTION: Full-length, untagged human recombinant histone H2A (residues 1-130; Genbank Accession #NM_021052.2; MW = 14.1 kDa) expressed in *E. coli*. Full-length, untagged human recombinant histone H2B (residues 1-126; Genbank Accession # NM_080593.2; MW = 13.9 kDa) expressed in *E. coli*.

PURITY: >95% by SDS-PAGE

SUPPLIED AS: _ µg/µL in 10 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM EDTA, 5 mM 2-mercaptoethanol, 10% glycerol

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (16% acrylamide) of 1µg and 2µg of RBC H2A-H2B dimer. Histones migrate anomalously at 17kDa (H2A) and 16kDa (H2B). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com