

## Product Information

<b>Description:</b>	Human recombinant histone octamer consisting of 2 molecules each of histones H2A encompassing amino acids 2-130(end), H2B encompassing amino acids (2-126(end), H3 encompassing 2-136(end), and H4 encompassing amino acids 2-103(end). Each recombinant histone has an N-terminal His-tag (6xHis). The recombinant complex was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	H2A (His-2-130(end)) / H2B (His-2-126(end)) / H3 (His-2-136(end)) / H4 (His-2-103(end))
<b>Concentration:</b>	3.34 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	45 mM Tris-HCl, pH 7.4, 2 M NaCl, and 10% glycerol
<b>MW:</b>	H2A: 15 kDa; H2B: 15 kDa; H3: 16 kDa; and H4: 12 kDa
<b>Genbank Accession:</b>	H2A: NM_033445; H2B: NM_003528; H3: NM_003532; H4: NM_003548
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

