

# Recombinant Nucleosome (*E. coli*-derived), Biotin-labeled, His-Tag

Catalog: 52048

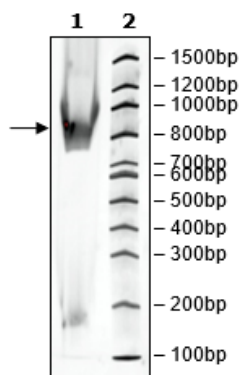
Lot: 220329-G

## Product Information

<b>Construct:</b>	Histone 2A (His-2-130(end)) / Histone H2B (His-2-126(end)) / Histone H3 (His-2-136(end)) / Histone H4 (His-2-103(end))-(Biotin)
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	0.90 mg/ml (by Bradford)
<b>Species:</b>	Human
<b>Formulated In:</b>	20 mM Tris-HCl pH 7.5, 1 mM DTT, 1 mM EDTA, and 20% glycerol
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	Histone H2A: NM_033445; Histone H2B: NM_003528; Histone H3: NM_003532; Histone H4: NM_003548
<b>MW:</b>	Histone H2A: 15 kDa; Histone H2B: 15 kDa; Histone H3: 16 kDa; Histone H4: 12 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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



### Biotin-Avidin Pulldown

