

## Data Sheet

### Histone Octamer, Biotin-labeled

Human, recombinant

**Catalog #:** 52061

**Lot #:** 131205-B      **Conc.:** 1.8 mg/ml

**Formulated in:** 45 mM Tris pH 7.5, 2 M NaCl, and 10% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Miller, K.M., Jackson, S.P., *Biochem. Soc. Trans.* 2012; **40(2)**: 370-6.
2. Raut, V.V., *et al. Ann. Bot.* 2011; **108(7)**:1235-46.

**Description:** Human recombinant histone octamer consisting of 2 molecules each of N-terminal His-tagged histones H2A (a.a. 2-130(end)), H2B (a.a. 2-126(end)), H3 (a.a. 2-137(end)), and H4 (a.a. 2-103(end)), expressed in *E. coli*. GenBank Accession Nos. are NM\_033445, NM\_003528, NM\_003532, and NM\_003548, respectively. H2B has been biotin-labeled via thiol modification of an additional C-terminal cysteine prior to assembly. Global MW = 117 kDa.

**Application:** A good substrate for histone methyltransferases.

## Quality Assurance

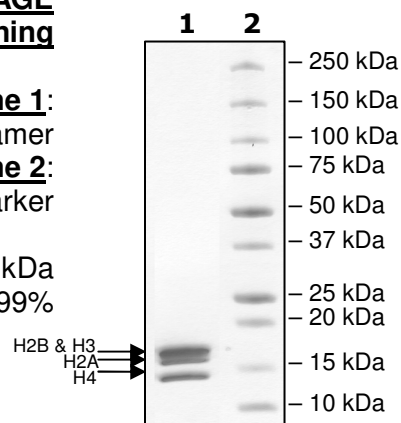
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
4  $\mu\text{g}$  Octamer

**Lane 2:**  
Protein Marker

**Global MW:** 117 kDa

**Purity:**  $\geq 99\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)