

## Data Sheet

### Histone H2A, His-Tag, Biotin-Labeled

Human, Recombinant, N-terminal His-tag,  
 Biotin Label at C-terminal Cysteine

**Catalog #:** 52049

**Lot#:** 130614      **Conc.:** 0.54 mg/ml

**Formulated in:** 8 mM PBS pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol and 3 mM DTT.

**Stability:** At least 6 months at -80°C  
 12 months from date of receipt, when stored as directed. Kit components require different storage conditions. Be sure to store each component at the proper temperature upon arrival.

**References:**

1. Osley, M.A. *Brief Funct. Genomic Proteomic* **5(3)**:179-89 (2006).
2. Wyrick, J.J. and Parra, M.A.. *Biochim Biophys Acta*. **1789(1)**:37-44 (2009).

**Description:**

Biotinylated Human Histone 2A, GenBank Accession No. NM\_033445, a.a. 2-130(end) with N-terminal His-tag and C-terminal Cys. MW = 15 kDa, expressed in an *E. coli* expression system. This protein was biotinylated using cysteine specific biotinylation at the C-terminus. Biotinylation was confirmed using western blot analysis with Strep-HRP.

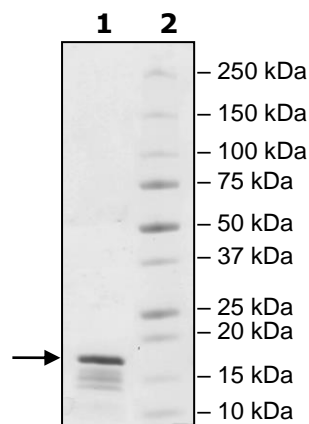
### **Quality Assurance**

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

**Lane 1:**  
 2 µg Histone  
 H2A+Cys

**Lane 2:**  
 Protein Marker

**MW:** 15 kDa  
**Purity:** ≥56%



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)



6044 Cornerstone Ct. West, Ste. E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MHHHHHSGRGKQGGKARAKAKSRSSRAGLQFPVGRVHRLLRKGNYSERVGAGAPV  
YLAHVLEYLTAEILELAGNAARDNKKTRIIPRHLQLAIRNDEELNKLLGRVTIAQGGVLPNI  
QAVLLPKKTESHHKAKGKC

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)