

Data Sheet

Histone H2B (full length)

Human, recombinant, N-terminal His-tag

Catalog #: 52026

Lot #: 110415 **Conc.:** 1.3 mg/ml

Formulated in: 50 mM Tris-HCl, pH 8.0, 150 mM NaCl, 10% glycerol, 1 mM PMSF and 1 mM DTT.

Stability: >6 months at -80°C

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).
3. Emre, N.C. and Berger, S.L. *Cold Spring Harb Symp Quant Biol*. **69**: 289-99 (2004).

Description:

Human Histone B, GenBank Accession No. NM_003528, a.a. 2-126(end) with N-terminal His-tag, MW = 14.6 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

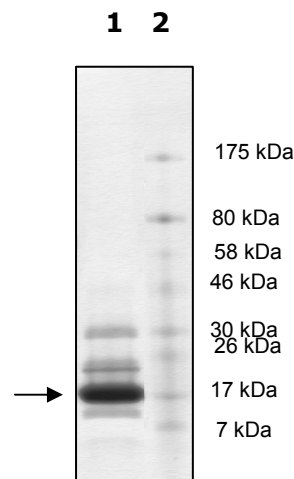
Quality Assurance

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

Lane 1:
Histone H2B

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 14.6 kDa
Purity: >65%



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

Please visit our website at: www.bpsbioscience.com