

Data Sheet

Histone H4 (full length)

Human, recombinant, N-terminal His-tag

Catalog #: 52028

Lot#: 130201 Conc.: 1.5 mg/ml

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

Stability: >6 months at -80°C

References:

1. Kornberg, R.D. (1977) *Annu. Rev. Biochem.*, **46**, 931-954.
2. van Holde, K.E. (1989) *Chromatin*, 1-497.

Description:

Human Histone 4 (HIST2H4A) (GenBank Accession No. NM_003548), a.a. 2-104(end) with N-terminal His-tag, MW = 12.1 kDa, expressed in an *E. coli* expression system.

Application:

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

Quality Assurance

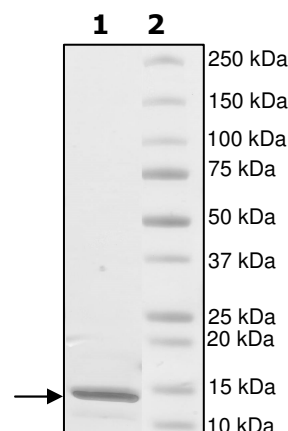
4-20 % SDS-PAGE Coomassie staining

Lane 1:
2 µg Histone H4

Lane 2:
Protein Marker

MW: 12.1 kDa

Purity: ≥93%



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