

Data Sheet

Histone H4 (2-58)

Human, recombinant, N-terminal His-tag

Catalog #: 52019

Lot#: 111201 Conc.: 0.68 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 16 mM
Glutathione, 20% Glycerol, and 3 mM DTT.

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 (GenBank Accession No. NM_003548), (2-58) with N-terminal GST-tag, MW = 32 kDa, expressed in an *E. coli* expression system.

Application:

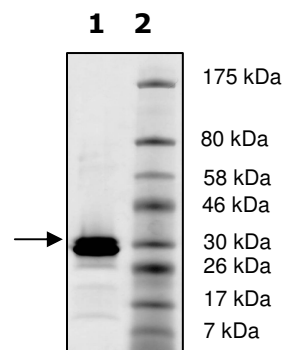
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2.4 µg Histone H4
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 32 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