

Data Sheet

Histone H3/H4 Tetramer

Human, recombinant, N-terminal His-tag

Catalog #: 52064

Lot#: 130926 **Conc.:** 0.82 mg/ml

Formulated in: 45 mM Tris pH 7.5, 2M NaCl, 1mM EDTA, 1mM DTT and 10% glycerol.

Stability: >6 months at -70°C

References:

1. Huang, C., *et al.*, *PLoS Genet.* 2013; **9(6)**: 1371.
2. Liu, W.H., *et al.*, *Nucleic Acids.* 2012; **40(22)**:11229-39.

Description:

Human recombinant histone tetramer, full length, consisting of 2 molecules each of histones H3 (a.a. 1-136(end)), and H4 (a.a. 1-103(end)), each with an N-terminal His-tag, expressed in an *E. coli* expression system. GenBank Accession Nos. are NM_003532, and NM_003548, respectively. Global MW = 57 kDa.

Application:

A perfect substrate for histone methyltransferases.

Quality Assurance

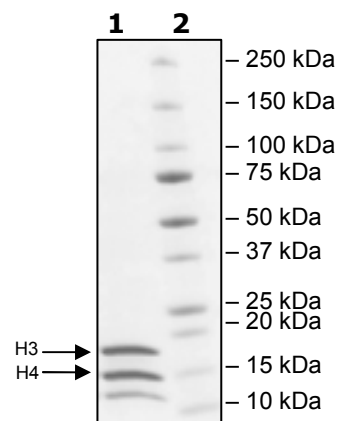
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg Tetramer

Lane 2:
Protein Marker

Global MW: 57 kDa

Purity: ≥80%



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequences:

Histone H3

MMHHHHHHHARTKQTARKSTGGKAPRKQLATKAARKSAPATGGVKKPHRYRPGTVALRE
IRRYQKSTELLIRKLPFQRLVREIAQDFKTDLRFQSSAVMALQEACEAYLVGLFEDTNLCAI
HAKRVTIMPKDIQLARRIRGERA

Histone H4

MGHHHHHHHSGRGKGGKGLGKGGAKRHRKVLRDNIQGITKPAIRRLARRGGVKRISGLIY
EETRGVLKVFLENVIRDAVITYTEHAKRKTVTAMDVVYALKRQGRTLYGFGG

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