

Data Sheet

Histone H1 (full length)

Human, recombinant, N-terminal His-tag

Catalog #: 52043

Lot#: 130521

Conc.: 1.57 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. Doenecke, D., *et al.*, *J. Mol. Biol.* 1986 Feb 5; **187(3)**:461-4.
2. Vyas, P., *et al.*, *J. Biol. Chem.* 2012 Apr 6; **287(15)**:11778-87.

Description:

Human Histone H1, also known as H1F0, (GenBank Accession No. NM_005318), a.a. 2-194(end) with N-terminal His-tag, MW = 21.7 kDa, expressed in an *E. coli* expression system.

Application:

Substrate for histone methyltransferase and acetyltransferase assays. Ideal for screening small molecular inhibitors of histone modifying enzymes for drug discovery and HTS applications.

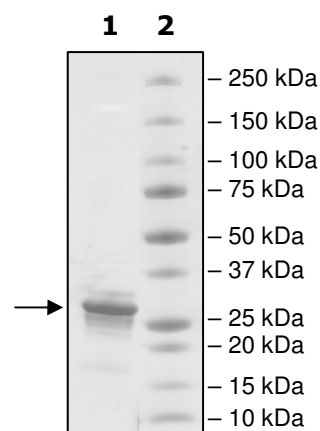
Quality Assurance

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

Lane 1:
2 µg Histone H1

Lane 2:
Protein Marker

MW: 22 kDa
Purity: ≥82%



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

Sequence:

MHHHHHHTENSTSAPAAKPKRAKASKKSTDHPKYSDMIVAAIQAEKNRAGSSRQSIQK
YIKSHYKVGENADSIKLSIKRLVTTGVLKQTKGVBGASGSFRLAKSDEPKKSVAFKTK
KEIKKVATPKKASKPKKAASKAPTCKKPKATPVKKAKKKLAATPKKAKKPKTVKAKPVK
ASKPKKAKPVKPKAKSSAKRAGKKK

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