

Data Sheet

Histone Octamer (full length)

Human, recombinant
Catalog #: 52040
Lot#: 130422 **Conc.:** 0.25 mg/ml

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

Stability: >6 months at -80°C

References:

1. Miller, K.M., Jackson, S.P., *Biochem Soc Trans.* 2012; **40(2)**: 370-6.
2. Raut, V.V., *et al.*, *Ann Bot.* 2011; **108(7)**:1235-46.

Description:

Human recombinant histone octamer consisting of 2 molecules each of histones H2A (a.a. 1-130(end)), H2B (a.a. 1-126(end)), H3 (a.a. 1-136(end)), and H4 (a.a. 1-103(end)), expressed in an *E. coli* expression system. GenBank Accession Nos. are NM_033445, NM_003528, NM_003532, and NM_003548, respectively. Global MW = 110 kDa.

Application:

A perfect substrate for histone methyltransferases.

Quality Assurance

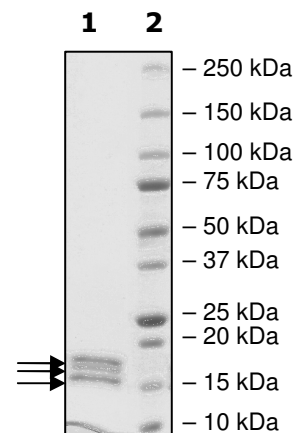
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 µg Octamer

Lane 2:
 Protein Marker

Global MW: 110 kDa
Purity: ≥99%

Histone H2B & Histone H3
 Histone H2A
 Histone H4



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 H2A

MSGRGKQGGKARAKAKSRSSRAGLQFPVGRVHRLLRKGNYSERVGAGAPVYLAADVLEY
LTAEILELAGNAARDNKKTRIIPRHLQLAIRNDEELNLLGRVTIAQGGVLPNIQAVLLPKKT
ESHHKAKGK

Histone H2B

MPEPAKSAPAPKKGSKKAVTKAQKKGKRRKRSRKESYSIYVYKVLKQVHPDTGISSKA
MGIMNSFVNDIFERIAGEASRLAHYNKRSTITSREIQTAVRLLLPGELAKHAVSEGTKAVTK
YTSSK

Histone H3

MARTKQTARKSTGGKAPRKQLATKAARKSAPATGGVKKPHRYRPGTVALREIRRYQKST
ELLIRKLPFQRLVREIAQDFKTDLRFQSSAVMALQEACEAYLVGLFEDTNLCAIHAKRVTIM
PKDIQLARRIRGERA

Histone H4

MGSGRGKGGKGLGKGGAKRHRKVLDRDNIQGITKPAIRRLARRGGVKRISGLIYEETRGVL
KVFLENVIRDAVITYTEHAKRKTVTAMDVVYALKRQGRTLYGFGG

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