

Data Sheet

Histone Octamer, H3K27C

Human, recombinant, N-terminal His-tag
Catalog #: 52088
Lot#: 140214 **Conc.:** 0.23 mg/ml

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

Stability: >6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Simon, M.D., *et al.* *Cell* 2007; **128**: 1003-1012.
2. Raut, V.V., Pandey, S.M., Sainis, J.K. *Ann Bot.* 2011; **108(7)**:1235-46.

Description:

Human recombinant histone tetramer 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. Each histone protein has an N-terminal His-tag. Histone H3 has K27C, C97A, and C111A mutations and was alkylated to produce the lysine analog, aminoethylcysteine. GenBank Accession Nos. are NM_033445, NM_003528, NM_003532, and NM_003548, respectively. Global MW = 116 kDa.

Application:

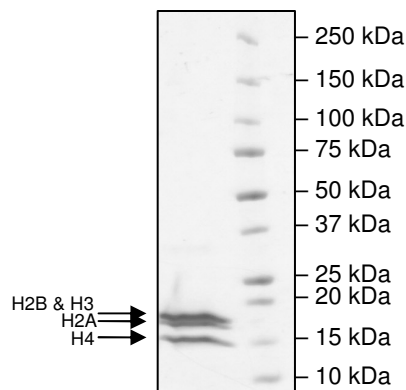
A perfect substrate for histone methyltransferases.

Quality Assurance

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

Lane 1:
4 μg Octamer
H3K27C-Unmodified
Lane 2:
Protein Marker

Global MW: 116 kDa
Purity: $\geq 90\%$



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