

Data Sheet

Histone H3-H4 Tetramer, H3K27C

Human, recombinant, N-terminal His-tag

Catalog #: 52084

Lot#: 140214 **Conc.:** 0.45 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 H3 (a.a. 2-136(end)) and H4 (a.a. 2-103(end)), expressed in an *E. coli* expression system. Histone H3 has K27C, C97A, and C111A mutations and was alkylated to produce the lysine analog, aminoethylcysteine. GenBank Accession Nos. NM_003532, and NM_003548, respectively. Each histone protein has an N-terminal His-tag. Global MW = 57 kDa.

Application:

A perfect substrate for histone methyltransferases.

Quality Assurance

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

Lane 1:

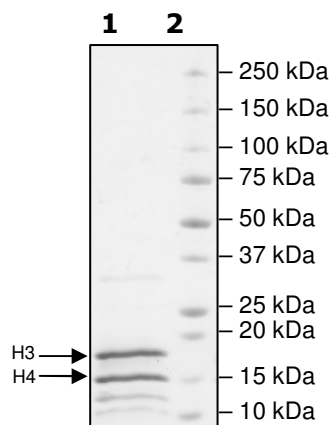
4 μg H3-H4 Tetramer
H3K27C-Unmodified

Lane 2:

Protein Marker

Global MW: 57 kDa

Purity: $\geq 80\%$



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