

## Data Sheet

### **Histone H3 K27C, Full Length**

Human, recombinant, N-terminal His-tag

**Catalog #:** 52080

**Lot#:** 140207

**Conc.:** 0.53 mg/ml

**Formulated in:** PBS (8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl), 20% glycerol.

**Stability:** >6 months at  $-80^{\circ}\text{C}$ . 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 H3 (a.a. 1-136(end)) with K27C, C97A, and C111A mutations, and an N-terminal His-tag, expressed in an *E. coli* expression system. GenBank Accession No. NM\_003532. The protein has been alkylated to produce the lysine analog, aminoethylcysteine. MW = 16 kDa.

**Application:**

A perfect substrate for histone methyltransferases.

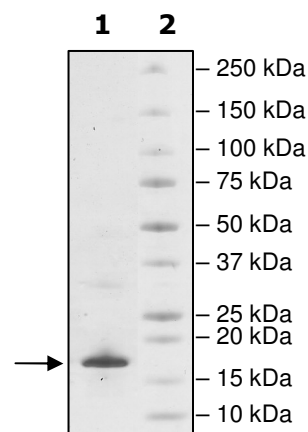
## **Quality Assurance**

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

**Lane 1:**  
2  $\mu\text{g}$  Histone H3K27C,  
Unmethylated

**Lane 2:**  
Protein Marker

**MW:** 16 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)