

## Data Sheet

### Histone Octamer, H3K27cMe3

Human, recombinant, N-terminal His-tag

**Catalog #:** 52091

**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  $-80^{\circ}\text{C}$ . 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 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. 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, trimethyl 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 and demethylases.

## Quality Assurance

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

**Lane 1:**

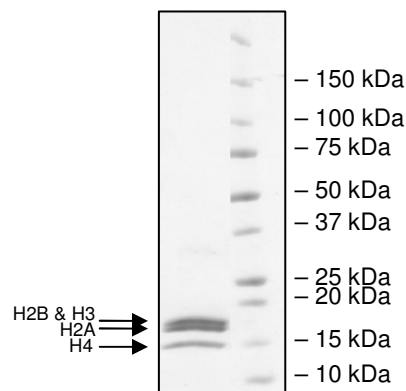
4  $\mu\text{g}$  Octamer  
 H3K27Cme3

**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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)