

## Data Sheet

### **JARID1C (KDM5C)**

Human recombinant, with N-terminal FLAG-tag and His-tag

**Catalog #:** 50112

**Lot #:** 140709      **Conc.** 1.3 mg/ml

**Formulated in:** 40 mM Tris-HCl, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 80 µg/mL and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

#### **References:**

1. Xu, J., *et al.* (2008). *PLoS One* **3(7)**:e2553.
2. Iwase, S., *et al.* (2007). *Cell* **128(6)**:1077-88.

**Description:** Human JARID1C, also known as SMCX (Smcy homolog, X-linked), KDM5C, or XE169, GenBank Accession No. NM\_004187), full length [a.a. 2-1560(end)] with N-terminal FLAG- and His-tags, MW= 181 kDa, expressed in Sf9 cells via a baculovirus expression system.

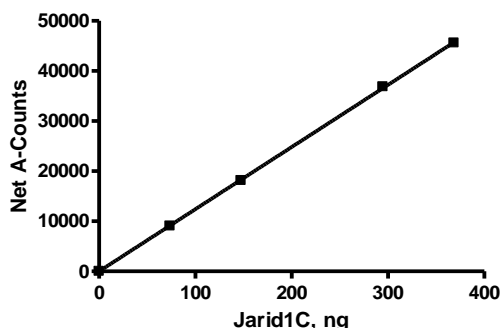
**Specific Activity:** 0.14 pmol/min/µg.

**Assay Conditions:** Assay was conducted using AlphaScreen® technology using BPS JARID1C Homogeneous Assay Kit, Catalog #50511. JARID1C enzyme is incubated with biotinylated histone H3 peptide substrate in JARID assay buffer in a 10 µl reaction for 2 hours at RT. AlphaLISA® anti-Mouse Acceptor beads (Perkin Elmer) are added, followed by the primary antibody addition and 30 min incubation at RT. AlphaScreen® Streptavidin-conjugated donor beads (Perkin Elmer) are added. After 10-15 minute incubation at room temperature, Alphacounts are measured.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

### **Specific Activity**



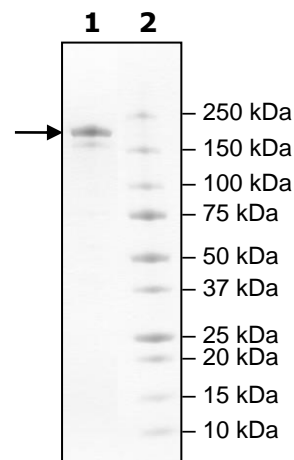
### **4-20% SDS-PAGE Coomassie staining**

**Lane 1:**  
2.2 µg JARID1C

**Lane 2:**  
Protein Marker

**MW:** 181 kDa

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

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