

## Data Sheet

### JARID1B (KDM5B), Full Length, FLAG-tag

Human, recombinant, N-terminal FLAG-Avi-tag

**Catalog #:** 50128

**Lot #:** 170511-3

**Conc.:** 0.18 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 80 µg/ml FLAG peptide, and 0.2 mM TCEP

**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:**

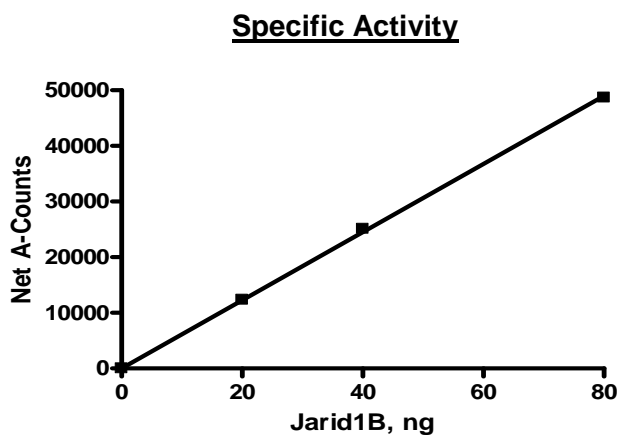
1. Yamane K, *et al.* (2007). *Mol. Cell* **25** (6): 801–12.
2. Barrett A, *et al.* (2007). *Int. J. Cancer* **121** (2): 265–75.

**Description:** Human JARID1B, also known as KDM5B and PLU-1 (GenBank Accession No. NM\_006618), a.a. 2-1544(end) with N-terminal FLAG-Avi-tag, MW=179 kDa, expressed in Sf9 cells via a baculovirus expression system.

**Specific Activity:** 0.11 pmol/min/µg  
 Assay conditions: 10 µl reaction mix containing 20 mM HEPES (pH 7.4), 50 mM NaCl, 500 µM  $\alpha$ -ketoglutarate, 25 µM iron, 2 mM ascorbic acid, 0.01% Tween-20, 1 mM TCEP, 0.5 µM biotinylated peptide substrate, and JARID1B. Incubate 1 hour at RT with slow shaking. Add antibody against demethylated product and anti-mouse acceptor beads. After 30 minutes, add Streptavidin-conjugated secondary antibody followed by Alpha Screening detection.

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

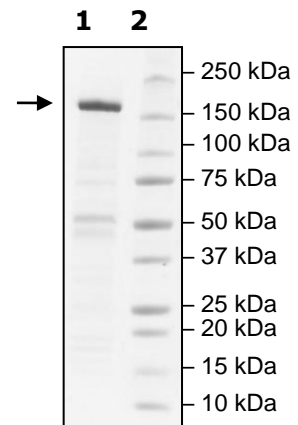
## Quality Assurance



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

**Lane 1:**  
 2.3 µg JARID1B  
**Lane 2:**  
 Protein Marker

**Purity:** 75%  
**MW:** 179 kDa



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)