

# JARID1B (KDM5B), Full Length, FLAG-Tag Recombinant

Catalog: 50128

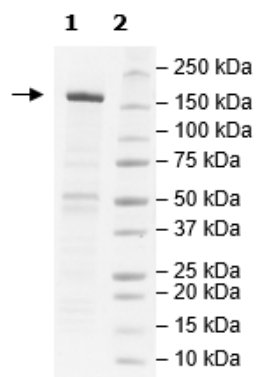
Lot: 150511-3

## Product Information

<b>Construct:</b>	JARID1B (KDM5B) (FLAG-Avi-2-1544(end))
<b>Concentration:</b>	0.18 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	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
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_006618
<b>MW:</b>	179 kDa
<b>Purity:</b>	75%
<b>Specific Activity:</b>	0.11 pmol/min/µg
<b>Assay Conditions:</b>	10 µl reaction mix containing assay buffer with 20 mM HEPES (pH 7.4), 50 mM NaCl, 500 µM α-ketoglutarate, 25 µM iron, 2 mM ascorbic acid, 0.01% Tween-20, 1 mM TCEP, 0.5 µM biotinylated peptide substrate, and JARID1 added to the wells. Incubate 1h at room temperature with slow shaking. Add antibody against demethylated K <sub>4</sub> peptide and anti-mouse acceptor beads. Incubate for 30 min, then add Streptavidin-conjugated secondary antibody followed by Alpha Screening detection.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



JARID1B Activity

