

JARID1B (KDM5B)

**CATALOG NO.:** PDM-21-360

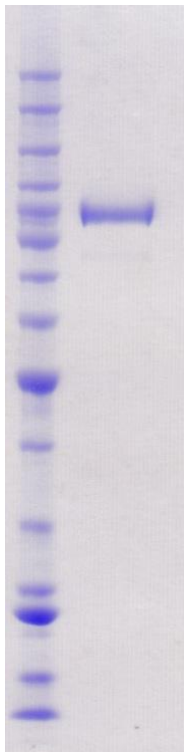
**LOT NO.:**

**DESCRIPTION:** Human recombinant JARID1B (residues 2-752; Genbank Accession # NM\_006618.3; MW = 90.7 kDa) expressed with an N-terminal FLAG tag in *Sf21* insect cells.

**PURITY:** >95% by SDS-PAGE

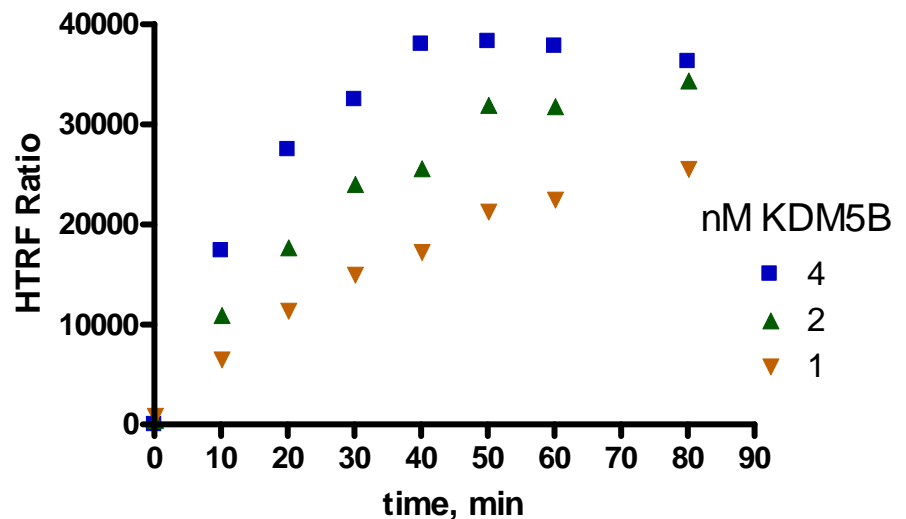
**SUPPLIED AS:**  $\mu\text{g}/\mu\text{L}$  in 50 mM Tris HCl, pH 7.5, 150 mM NaCl, 10% glycerol, 1 mM TCEP

**STORAGE:**  $-70^{\circ}\text{C}$ . Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4  $\mu\text{g}$  of RBC JARID1B (MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.

**Demethylase activity of KDM5B**



**KDM5B Demethylase Activity Assay.** B-H3(1-21)K4Me3 peptide demethylation was monitored using Cisbio HTRF detection (cat # 61KA2KAD and 610SAXLA). The 10  $\mu\text{L}$  reaction contained 20  $\mu\text{M}$  Fe(II), 20  $\mu\text{M}$   $\alpha$ -KG, 100  $\mu\text{M}$  ascorbate, 30 nM peptide and variable concentration of KDM5B. Reactions were quenched at specified time by the addition of detection reagent. Fluorescence emission (665 and 620 nm) was read using a Synergy H4 plate reader (Biotek) following 2h incubation.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.