

## Recombinant JMJD2D / KDM4D protein

---

**Catalog No:** 31459, 31859

**Expressed In:** Baculovirus

**Quantity:** 20 µg

**Concentration:** 0.25 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant JMJD2D / KDM4D protein is supplied at a concentration of 0.25 µg/µl in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** KDM4D (lysine (K)-specific demethylase 4D), also known as JMJD2D (Jumonji Domain Containing 2D) is a protein that functions as a histone demethylase that preferentially demethylates di- and trimethylated lysine 9 residues of histone H3, while it has no activity on monomethylated H3K9 residues.

**Protein Details:** Recombinant JMJD2D / KDM4D (accession number NP\_060509.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 63.2 kDa. The recombinant protein is >60% pure by SDS-PAGE.

**Application Notes:** Recombinant JMJD2D / KDM4D is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

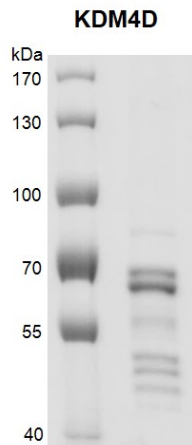
**Specific Activity:** H3K9me3 demethylase.

**Histone Demethylase Assay Conditions:**

50 mM HEPES pH 7.5, 0.02% Triton X-100, 100 µM 2OG, 100 µM Ascorbate, 50 µM (NH<sub>4</sub>)<sub>2</sub>Fe(SO<sub>4</sub>)<sub>2</sub>·6H<sub>2</sub>O, 1 mM TCEP, 100 nM Recombinant JMJD2D / KDM4D protein, and 3.3 µM H3K9me3 (aa 1-21) peptide at 2 hours at room temperature. MALDI-TOF was used for detection.

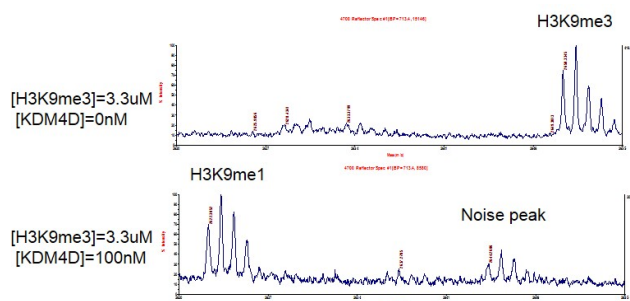
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



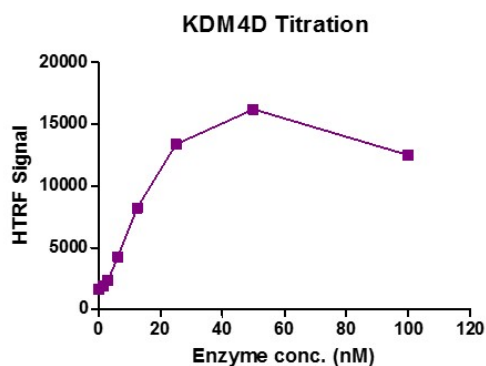
#### Recombinant JMJD2D / KDM4D protein gel.

JMJD2D / KDM4D protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.



#### MALDI-TOF Assay for JMJD2D / KDM4D activity .

3.3  $\mu$ M H3K9me3 peptide was incubated with 100 nM JMJD2D / KDM4D in reaction buffer for 2 hours at room temperature. The reaction product was detected by MALDI-TOF. Single 3.3  $\mu$ M H3K9me3 peptide was used as a negative control. Catalytic Ability: >66 turnovers / enzyme molecule.



#### HTRF Assay for Recombinant JMJD2D / KDM4D activity.

3.3  $\mu$ M H3K9me3 peptide was incubated with JMJD2D / KDM4D protein in reaction buffer for 1 hour at room temperature. Anti-H3K9me2 antibody was used to detect reaction product.