

## Recombinant FBXL10 / KDM2B protein

---

**Catalog No:** 31455, 31855

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant FBXL10 / KDM2B protein is supplied in 25 mM HEPES-NaOH, pH 7.5, 300 mM NaCl, 20% glycerol, 0.04 % Triton X-100, 0.5 mM TCEP.

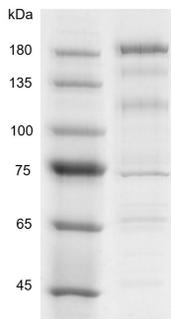
**Background:** **KDM2B (lysine (K)-specific demethylase 2B)**, also known as **FBXL10 (F-box and leucine-rich repeat protein 10)** is a histone demethylase that preferentially demethylates trimethylated lysine 4 (K4me3) and dimethylated lysine 36 (K36me2) of histone H3. KDM2B displays weak or no activity for mono- and trimethylated H3K36. KDM2B preferentially binds the transcribed region of ribosomal RNA and represses the transcription of ribosomal RNA genes which results in inhibition of cell growth and proliferation. KDM2B may also serve as a substrate-recognition component of the SCF (SKP1-CUL1-F-box protein)-type E3 ubiquitin ligase complex.

**Protein Details:** Recombinant FBXL10 / KDM2B was expressed in a baculovirus expression system as the full length proteins (accession number NP\_115979.3) with an N-terminal 6xHis tag and FLAG-Tag. The molecular weight of FBXL10 / KDM2B is 158.2 kDa. The recombinant protein is >60% pure by SDS-PAGE.

**Application Notes:** This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

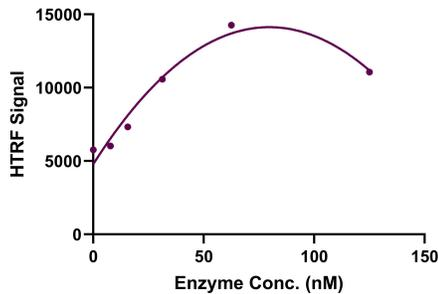
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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

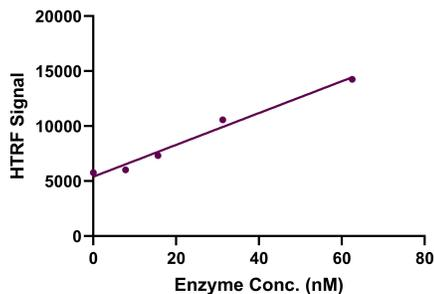
**FBXL10 / KDM2B**

**Recombinant FBXL10 / KDM2B protein**

7.5% SDS-PAGE Coomassie staining

MW: 158.2 kDa

 Purity:  $\geq 60\%$ 
**KDM2B Titration**

**HTRF assay for FBXL10 / KDM2B activity**

1  $\mu\text{M}$  H3K36me2 peptide was incubated with different concentrations of FBXL10 / KDM2B protein in 10  $\mu\text{l}$  reaction system containing 50 mM HEPES pH 7.5, 1 mM TCEP, 0.02% Triton X100, 100  $\mu\text{M}$  2-oxoglutarate, 100  $\mu\text{M}$  ascorbate, 50  $\mu\text{M}$   $(\text{NH}_4)_2\text{Fe}(\text{SO}_4)_2 \cdot 6\text{H}_2\text{O}$  and for 1 hours at room temperature then, then 10  $\mu\text{l}$  antiH3K36me1 antibody and SA-XL665 mixture (1:100 dilution in the same buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

**KDM2B Titration**

**HTRF assay for FBXL10 / KDM2B activity**

1  $\mu\text{M}$  H3K36me2 peptide was incubated with different concentrations of FBXL10 / KDM2B protein in 10  $\mu\text{l}$  reaction system containing 50 mM HEPES pH 7.5, 1 mM TCEP, 0.02% Triton X100, 100  $\mu\text{M}$  2-oxoglutarate, 100  $\mu\text{M}$  ascorbate, 50  $\mu\text{M}$   $(\text{NH}_4)_2\text{Fe}(\text{SO}_4)_2 \cdot 6\text{H}_2\text{O}$  and for 1 hours at room temperature then, then 10  $\mu\text{l}$  antiH3K36me1 antibody and SA-XL665 mixture (1:100 dilution in the same buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.