

## Recombinant KDM2A / FBXL11 protein

---

**Catalog No:** 31485, 31885

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant FBXL11 / KDM2A protein was expressed in Sf9 cells and is supplied in 25 mM HEPES pH 7.5, 500 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

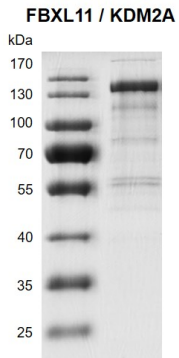
**Background:** **Lysine (K)-specific demethylase 2A (KDM2A)**, also known as FBXL11, JHDM1A, CXXC8, nucleates at CpG islands and specifically demethylates 'Lys-36' of histone H3, thereby playing a central role in histone code. Preferentially demethylates dimethylated H3 'Lys-36' residue while it has weak or no activity for mono- and trimethylated H3 'Lys-36'. May also recognize and bind to some phosphorylated proteins and promote their ubiquitination and degradation. Required to maintain the heterochromatic state. Associates with centromeres and represses transcription of small non-coding RNAs that are encoded by the clusters of satellite repeats at the centromere. Required to sustain centromeric integrity and genomic stability, particularly during mitosis.

**Protein Details:** Recombinant KDM2A / FBXL11 (accession number NP\_036440.1) was expressed in Sf9 and contains an N-terminal FLAG-Tag with a molecular weight of 134 kDa. The recombinant protein is >75% pure by SDS-PAGE.

**Application Notes:** Recombinant KDM2A / FBXL11 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling. Recombinant KDM2A / FBXL11 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**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 KDM2A / FBXL11 protein gel.**

KDM2A / FBXL11 protein was run on a 8% SDS-PAGE gel and stained with Coomassie blue.

MW: 128 kDa

Purity: > 85%

**MALDI-TOF for KDM2A / FBXL11 protein activity**

3  $\mu$ M H3K36me2 (21-44 aa) peptide was incubated with 0, 100 nM of KDM2A / FBXL11 protein in reaction buffer containing 50 mM HEPES-NaOH pH 7.5, 100  $\mu$ M 2-oxoglutarate, 100  $\mu$ M ascorbate, 50  $\mu$ M (NH<sub>4</sub>)<sub>2</sub>Fe(SO<sub>4</sub>)<sub>2</sub>·6H<sub>2</sub>O, 1 mM TCEP for 2 hr at room temperature. Single H3K36me2 peptide was used as negative control. MALDI-TOF was used for detection.

