

## Recombinant KDM1B / LSD2 protein



**Catalog No:** 31479, 31879

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant KDM1B / LSD2 protein is supplied in 25 mM Tris, pH 8, 300 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

**Background: KDM1B / LSD2** (Lysine (K)-Specific Demethylase 1B / Lysine-Specific Histone Demethylase 2) is a histone demethylase that demethylates 'Lys-4' of histone H3, a specific tag for epigenetic transcriptional activation, thereby acting as a corepressor. Required for *de novo* DNA methylation of a subset of imprinted genes during oogenesis. Demethylates both mono- and di-methylated 'Lys-4' of histone H3. Has no effect on tri-methylated 'Lys-4', mono-, di- or tri-methylated 'Lys-9', mono-, di- or tri-methylated 'Lys-27', mono-, di- or tri-methylated 'Lys-36' of histone H3, or on mono-, di- or tri-methylated 'Lys-20' of histone H4 (by similarity).

**Protein Details:** Full-length Recombinant KDM1B / LSD2 (accession number XM\_005248926.1) was expressed in Sf9 cells and contains an N-terminal GST-Tag with an observed molecular weight of 116 kDa. The recombinant protein is >92% 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.

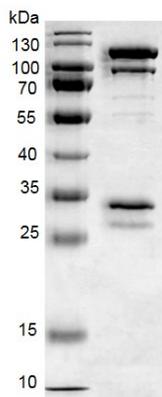
### References:

This product was used in the following publications:

*J. Biol. Chem.* (2016). 291(47): 24756-24767. PMID: 27634040.

**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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

**LSD2 / KDM1B**



**Recombinant KDM1B / LSD2 protein gel.**

Recombinant KDM1B / LSD2 run on an 8% SDS-PAGE gel and stained with Coomassie blue.