

## KDM1A Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # K421-31G**

Lot # 1158-2

### Product Description

Recombinant human KDM1A (172-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC048134](#).

### Gene Aliases

AOF2; BHC110; LSD1; KIAA0601

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

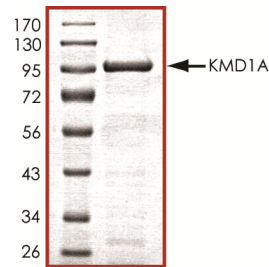
### Scientific Background

KDM1A or lysine (K)-specific demethylase 1A is a nuclear protein containing a SWIRM domain, a FAD-binding motif, and an amine oxidase domain that is a component of several histone deacetylase complexes and act as a histone demethylase. (1) H3K4 histone demethylase activity of KDM1A is partly responsible for the repressive activity of TAL1 and restricts TAL1 function in hematopoiesis (1). KDM1A plays an essential role for CoREST in demethylation of H3K4 both in vitro and in vivo (2).

### References

- Hu, X. et.al : LSD1-mediated epigenetic modification is required for TAL1 function and hematopoiesis. Proc. Nat. Acad. Sci. 106: 10141-10146, 2009.
- Lee, M. G. et.al : An essential role for CoREST in nucleosomal histone 3 lysine 4 demethylation. Nature 437: 432-435, 2005

### Purity



The purity of KDM1A was determined to be **>95%** by densitometry. Approx. MW **100kDa**.

## KDM1A Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	K421-31G
Specific Lot Number	1158-2
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**