

KDM2A Protein

Recombinant human protein expressed in Sf9 cells

Catalog # K422-31G

Lot # J665-3

Product Description

Recombinant human KDM2A (1-748) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_012308](#).

Gene Aliases

CXXC8; FBL11; FBL7; FBXL11; JHDM1A; LILINA

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

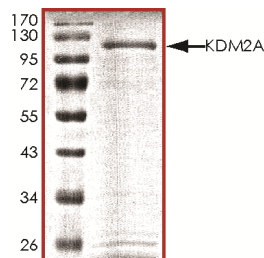
Scientific Background

KDM2A or lysine (K)-specific demethylase 2A is a member of the F-box protein family that are components of the modular E3 ubiquitin protein ligases called SCFs (SKP1-cullin-F-box), which function in phosphorylation-dependent ubiquitination (1). KDM2A belongs to the Fbls class and, in addition to the presence of an F-box, contains at least six highly degenerated leucine-rich repeats which play a role in epigenetic silencing. The demethylase activity of the JmjC domain-containing proteins is conserved from yeast to human (2).

References

1. Cenciarelli, C. et al: Identification of a family of human F-box proteins. *Curr. Biol.* 9: 1177-1179, 1999.
2. Tsukada, Y. et al: Histone demethylation by a family of JmjC domain-containing proteins. *Nature* 439: 811-816, 2006.

Purity



The purity of KDM2A was determined to be **>90%** by densitometry. Approx. MW **115 kDa**.

KDM2A Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number **K422-31G**

Specific Lot Number **J665-3**

Purity **>90%**

Concentration **0.1 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
www.signalchem.com

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