

KDM2B Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog # K422-30BG

Lot # P1809-6

Product Description

Recombinant full-length human KDM2B was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC115380](#).

Gene Aliases

FBXL10, CXXC2, Fbl10, JHDM1B, PCCX2

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.

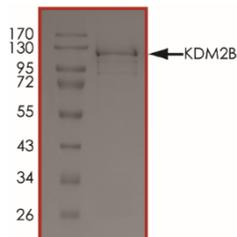
Scientific Background

KDM2B is a member of the F-box protein family which is characterized by an approximately 40 amino acid motif, the F-box proteins which constitute one of the four subunits of ubiquitin protein ligase complex called SCFs (SKP1-cullin-F-box), which function in phosphorylation-dependent ubiquitination (1). KDM2B functions in a DNA repair process other than mismatch repair. KDM2B is highly expressed in aggressive brain tumors, which play a main role in cancer development (2). KDM2B is a key regulator of JUN function.

References

1. Frescas, D. et al: JHDM1B/FBXL10 is a nucleolar protein that represses transcription of ribosomal RNA genes. *Nature* 450: 309-313, 2007.
2. Jin, J. et al: Systematic analysis and nomenclature of mammalian F-box proteins. *Genes Dev.* 18: 2573-2580, 2004.

Purity



The purity of KDM2B was determined to be **>75%** by densitometry. Approx. MW ~**120kDa**.

KDM2B Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	K422-30BG
Lot #	P1809-6
Purity	>75%
Concentration	0.05 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.