

L3MBTL1 3xMBT Domain, Recombinant Human

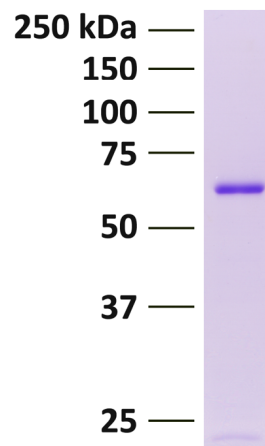
Catalog No. 15-0043
Lot No. 13289001
Pack Size 100 µg



Type MBT **Expressed In** *E. coli*
Mol. Wgt. 68 kDa **Epitope Tag** GST

Product Description:

Recombinant human L3MBTL1 (LMBL1, KIAA0681, L3MBT, L3MBTL, accession Q9Y468, amino acids 190 to 530) containing an N-terminal GST tag, expressed in *E. coli*. L3MBTL1 3xMBT contains the three MBT repeat domains present in the protein. L3MBTL1 is a member of the Polycomb group of proteins and binds to broad range of proteins that exhibit mono- and dimethylation, including histone H1, histone H4, p53 and Rb. The truncated version of L3MBTL1 present in the 3xMBT domain protein serves as a useful tool for the detection of mono- and dimethylation of a wide variety of proteins.



Protein Gel Data: L3MBTL1 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

L3MBTL1 3xMBT Domain, Recombinant Human is useful for the detection of protein mono- and dimethylation in far-Western assays, for the immunoprecipitation of methylated proteins in pull-down experiments, or as a readout of methyltransferase enzyme activity.

References:

Moore KE et al (2013) *Mol Cell* 50: 444-456.

This product is for *in vitro* research use only and is not intended for use in humans or animals.