

# L3MBTL1 3xMBT Domain, D355N, Recombinant Human

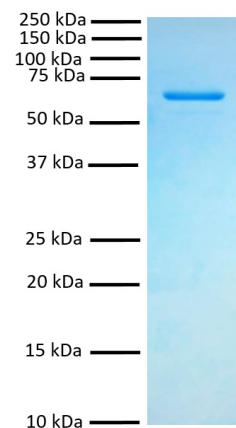
**Catalog No.** 15-0056  
**Lot No.** 16273001  
**Pack Size** 100 µg



**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 the wild type protein binds to broad range of proteins that exhibit mono- and dimethylation, including histone H1, histone H4, p53 and Rb. The D355N mutant version contains an asparagine at position 355 which abrogates methyl-lysine binding.



**Protein Gel Data:** L3MBTL1 3xMBT Domain, D355N, Recombinant Human (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 50 mM Tris-HCl pH 7.5, 300 mM NaCl, 1.5 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, D355N, Recombinant Human is useful as a non-binding control protein when studying wild type L3MBTL1.

## References:

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