

**L3MBTL4 (GST)**

**CATALOG NO.:** RD-11-404

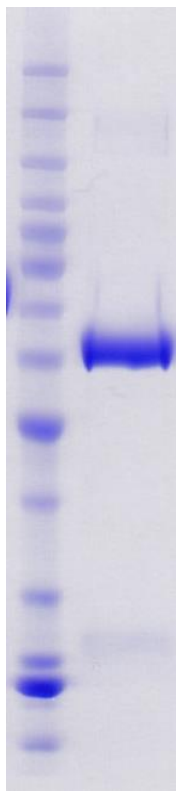
**LOT NO.:**

**DESCRIPTION:** Human recombinant L3MBTL4 (residues 49-376; Genbank Accession # NM\_173464.3; MW = 66.5 kDa) expressed as an N-terminal GST-fusion and C-terminal His-tag protein in *E. coli*.

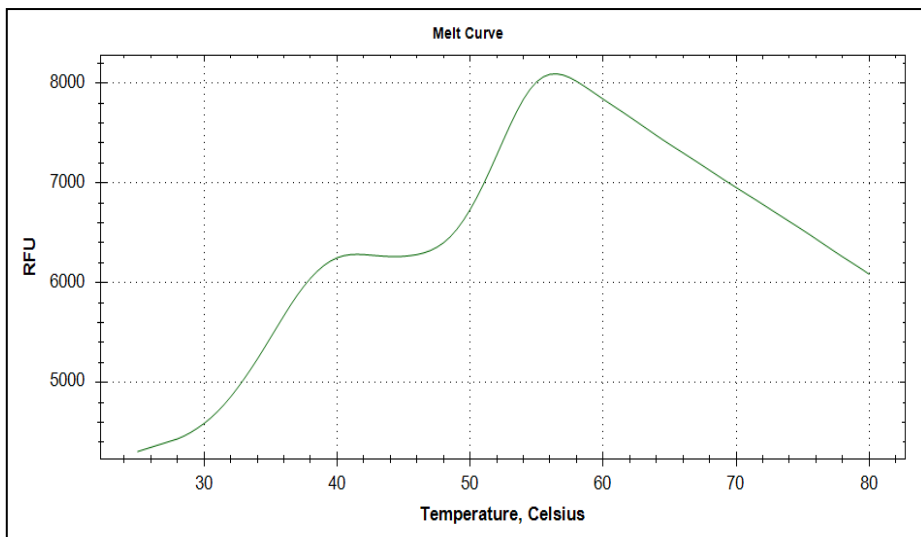
**PURITY:** >85% by SDS-PAGE

**SUPPLIED AS:**  $\_ \mu\text{g}/\mu\text{L}$  in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4  $\mu\text{g}$  of RBC L3MBTL4 (GST).** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Differential Scanning Fluorimetry of RBC L3MBTL4 (GST).** Thermal denaturation of L3MBTL4 (GST) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). The apo form of L3MBTL4 (GST) displays a Tm of 35.0°C.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355  
 requests@reactionbiology.com www.reactionbiology.com