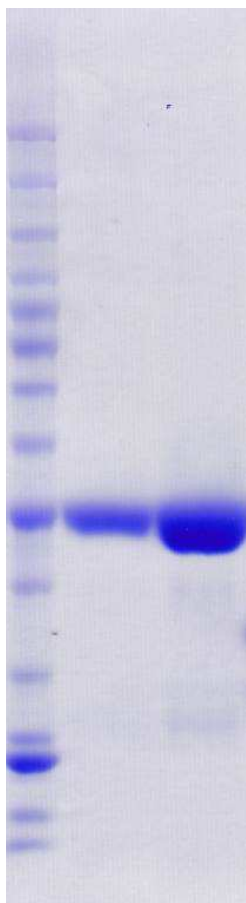


**SCML2 (GST)**
**CATALOG NO.:** RD-11-408

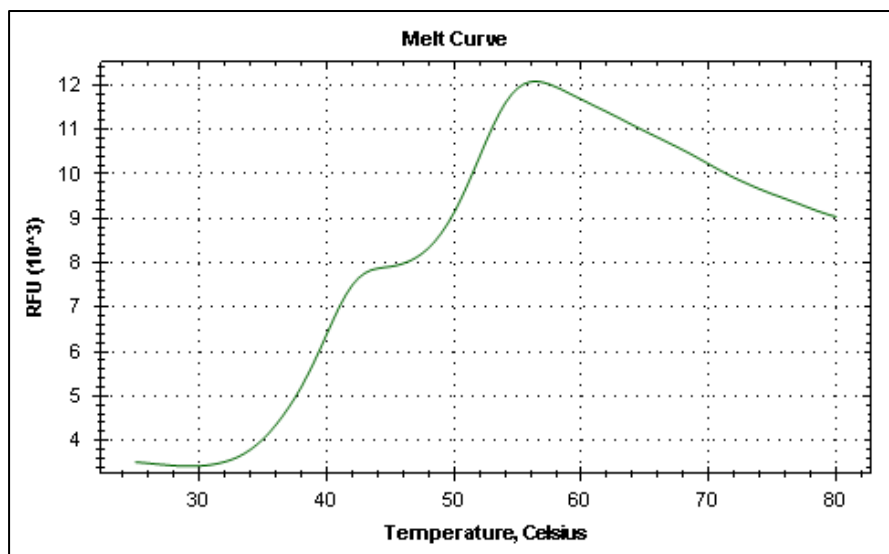
**LOT NO.:**
**DESCRIPTION:** Human recombinant SCML2 (residues 29-243; Genbank Accession # NM\_006089.2; MW = 51.2 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**PURITY:** >95% 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 as determined by OD<sub>280</sub>.

**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 10, 4  $\mu\text{g}$  of RBC SCML2 (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 SCML2 (GST).** Thermal denaturation of SCML2 (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 SCML2 (GST) displays a T<sub>m</sub> of 40.0°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