

Mras (wildtype) (GST)

CATALOG NO.: MSC-11-630

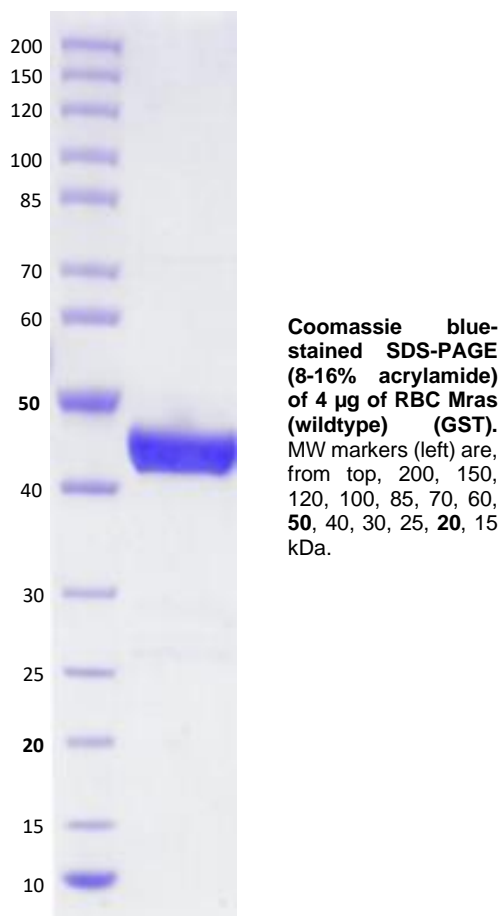
LOT NO.:

**DESCRIPTION:** Human recombinant Mras (wildtype) (residues 2-182; Genbank Accession #NM\_001085049.3 MW = 47.6 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**PURITY:** > 95% by SDS-PAGE.

**SUPPLIED AS:** \_\_\_ µg/µL in 25mM Tris-HCl pH 8.0, 500mM NaCl, 5% glycerol, 5mM MgCl<sub>2</sub>, 2mM DTT as determined by OD<sub>595</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 µl) or storage of diluted enzyme is not recommended.



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