

MIRO1 (Mitochondrial Rho GTPase 1)

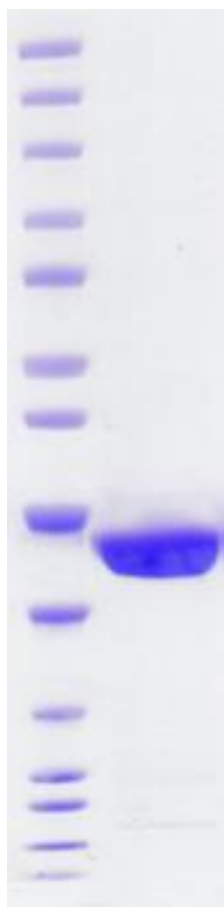
CATALOG NO.: MSC-11-682

DESCRIPTION: Human recombinant MIRO1 protein (residues 186-591; Genbank Accession # NM_018307.5; MW = 50.2 kDa) expressed as an N-terminal 8xHis-tag protein in *E.coli*.

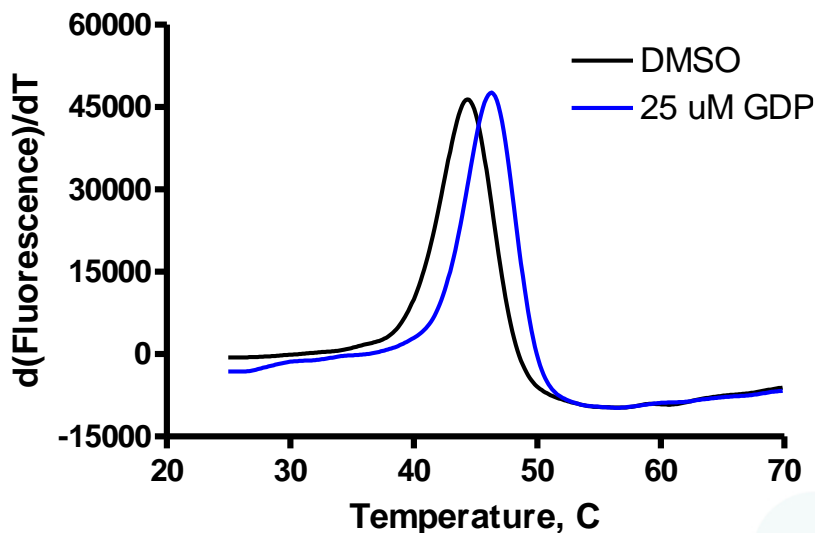
PURITY: > 95% by SDS-PAGE.

SUPPLIED AS: ___ $\mu\text{g}/\mu\text{L}$ in 25mM Tris-HCl pH 8, 500mM NaCl, 5mM MgCl₂, 5% glycerol, 0.25mM TCEP as determined by OD₅₉₅.

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.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC MIRO1. MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, **50**, 40, 30, 25, 20, 15, & 10 kDa.



Differential Scanning Fluorimetry of RBC MIRO1 in Presence or Absence of GDP. Thermal denaturation of MIRO1 is monitored on the QS5 384 QPCR SYSTEM (Thermo Fisher) by increased binding and fluorescence of GloMelt Thermal Shift dye (Biotium). Addition of 25 μM GDP stabilizes the protein and shifts the T_m (inflection point) from 44.3°C to 46.3°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