

MPP8-[CHR] (His) {M-phase phosphoprotein 8-[CHR] (His)}

CATALOG NO.: RD-11-462

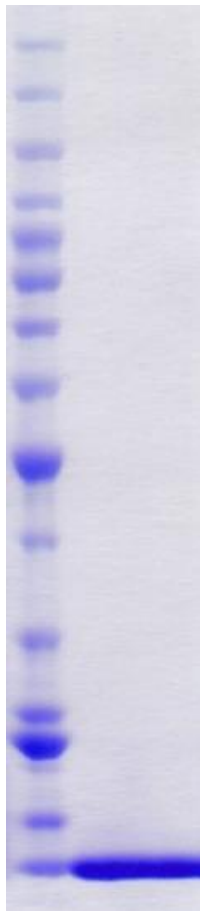
LOT NO.:

DESCRIPTION: Human recombinant MPP8-[CHR] (residues 55-116; Genbank Accession # NM_017520.3; MW = 9.9 kDa) expressed as an N-terminal His-tag fusion protein in *E. coli*.

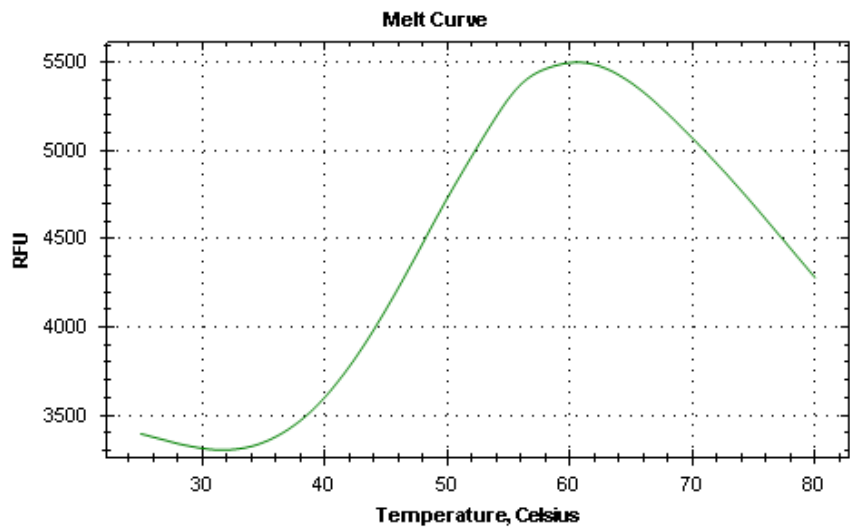
PURITY: >95% by SDS-PAGE

SUPPLIED AS: ___ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol as determined by OD₂₈₀.

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 MPP8-[CHR] (His). 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 MPP8-[CHR] (His). Thermal denaturation of MPP8-[CHR] (His) 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 MPP8-[CHR] (His) displays a T_m of 48.0°C.

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