

WRN (aa 500-946)

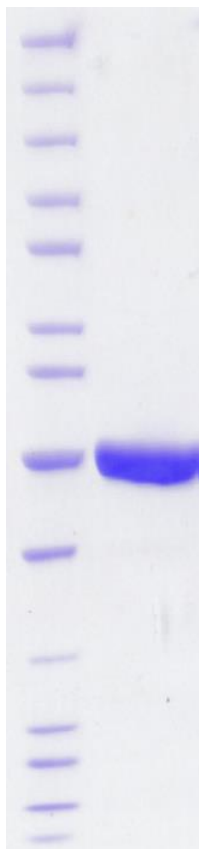
CATALOG NO.: HLC-21-674

DESCRIPTION: Human recombinant WRN protein (residues 500-946; Genbank Accession # AAF06162.1; MW = 52.5 kDa) expressed as a C-terminal His-tag protein in *SF9* cells.

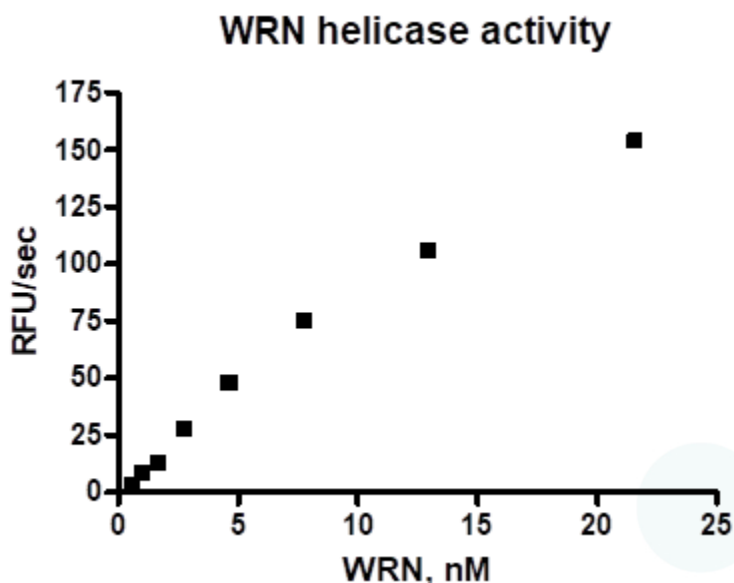
PURITY: > 95% by SDS-PAGE.

SUPPLIED AS: µg/µL in 50mM Tris-HCl, pH 7.5, 500mM NaCl, 10% glycerol, 0.25mM TCEP 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 of RBC WRN (aa 500-946). MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, 50, 40, 30, 25, 20, 15, & 10 kDa.



Helicase activity of WRN was assessed using a fluorescent forked DNA substrate (TAMRA/BHQ2 labels) in the presence of 4 mM ATP.

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