

**PIF1 (ATP-dependent DNA helicase PIF1)****CATALOG NO.:** HLC-11-679**DESCRIPTION:** Human recombinant PIF1 protein (residues 202-620; Genbank Accession # NM\_001286496.2; MW = 47.8 kDa) expressed with a C-terminal StrepII-tag protein in *E.coli*.**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 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 µl) or storage of diluted enzyme is not recommended.

**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC PIF1.** MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, **50**, 40, 30, 25, **20**, 15, & 10 kDa.

---

*Enzymatic activity has not been tested.*

---

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