

**BLM (RecQ-like DNA helicase BLM)**

**CATALOG NO.:** HLC-11-678

**DESCRIPTION:** Human recombinant BLM protein (residues 636-1298; Genbank Accession # NM\_000057.4; MW = 77.7 kDa) expressed with a C-terminal StrepII tag 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 BLM.** 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