



Product #MV100-10

**Purified Human Mitochondrial
RNA Polymerase**

h-mtRNAP; POLRMT

10 µg in 20 µl

Lot # 141130MV

Expiration:

Product Description: Human mitochondrial RNA polymerase (h-mtRNAP or **POLRMT**) is a 134 kDa, single-subunit DNA-dependent RNA polymerase resident in human mitochondria. POLRMT transcribes the mitochondrial genome to produce the mRNA's and tRNA's that drive mitochondrial protein synthesis, as well as RNA primers required for mitochondrial DNA replication.

Concentration: 0.5 mg/mL in 10 mM Tris (pH 8.0), 1 mM DTT, 100 mM NaCl and 20% glycerol.

Storage: Stable for 6 months at -80°C from date of shipment. Aliquot to avoid repeated freezing and thawing. Working stocks may be stored at -20°C.

Source & Purity: Recombinant POLRMT is purified from *E. coli* and does not incorporate an affinity tag. POLRMT shows apparent homogeneity by SDS-PAGE (500 ng of protein analyzed by 15% SDS- polyacrylamide gel electrophoresis and Coomassie blue staining).

