



Product #MV300-10

**Purified Human Mitochondrial
Transcription Factor B**

(h-mtTFB2)

10 µg in 20 µl

Lot # 141130MV3

Expiration:

Product Description: Human mitochondrial Transcription Factor B2 (h-mtTFB2) is a 42 kDa, single-subunit protein resident in human mitochondria that is needed for *in vivo* mitochondrial transcription. h-mtTFB2 can stimulate promoter specific mtDNA transcription in an *in vitro* assay system comprising purified mitochondrial RNA polymerase (h-mtRNAP) and mitochondrial transcription factor A (h-mtTFA). h-mtTFB2 is homologous to bacterial rRNA methyltransferases.

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: h-mtTFB2 is a recombinant purified from *E. coli*; it does not incorporate an affinity tag. h-mtTFB2 shows apparent homogeneity by SDS-PAGE (500 ng of protein analyzed by 15% SDS- polyacrylamide gel electrophoresis and Coomassie blue staining).

