



Product #MV200-10

**Purified Human Mitochondrial
Transcription Factor A
(h-mtTFA; aka TFAM)**

10 µg in 20 µl

Lot # 141130MV2

Expiration:

Product Description: Human mitochondrial Transcription Factor A (h-mtTFA) is a 25 kDa, single-subunit protein resident in human mitochondria. It is a HMG-box DNA-binding protein that is essential for initiation of mitochondrial transcription from both the human mitochondrial light-strand promoter (LSP) and heavy strand promoter (HSP1).

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

