

TUFM Polyclonal Antibody

(Catalog # A52852)

Background

This protein promotes the GTP-dependent binding of aminoacyl-tRNA to the A-site of ribosomes during protein biosynthesis.

Description

TUFM Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P49411

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Elongation factor Tu, mitochondrial protein (46-290AA)

Storage

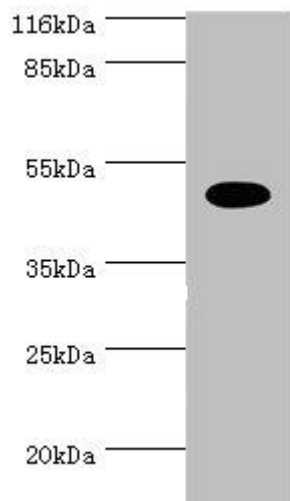
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

EF-Tu, P43, TUFM

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: TUFM Polyclonal Antibody at 2ug/ml+U251 whole cell lysate

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 50kDa

Observed band size: 50kDa