

MTMR7 Polyclonal Antibody

(Catalog # A56531)

Background

Phosphatase that acts on lipids with a phosphoinositol headgroup.

Description

MTMR7 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9Y216

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Myotubularin-related protein 7 protein (1-660AA)

Storage

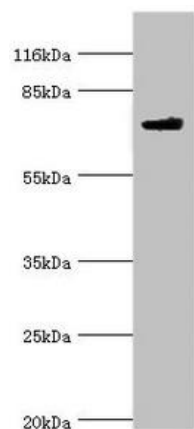
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

MTMR7

Application

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



Western blot

All lanes: MTMR7 antibody at 2ug/ml + Mouse brain tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 76, 65 kDa

Observed band size: 76 kDa