

RMDN1 Polyclonal Antibody

(Catalog # A66914)

Description

RMDN1 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

Q96DB5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Regulator of microtubule dynamics protein 1 protein (13-314AA)

Storage

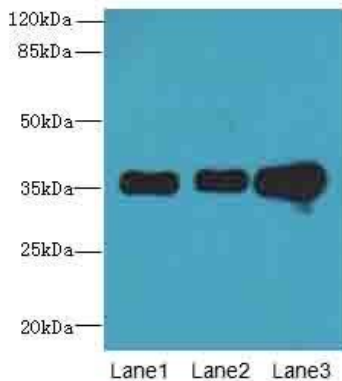
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Regulator of microtubule dynamics protein 1, RMD-1, hRMD-1, Protein FAM82B, RMDN1, FAM82B, CGI-90

Application

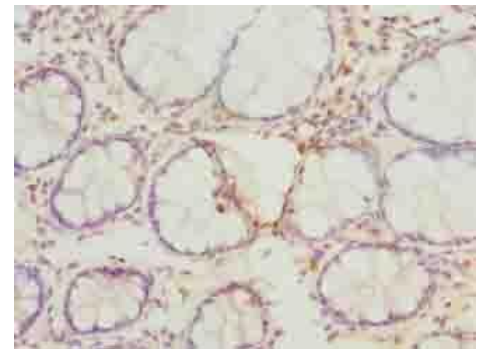
ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



WB: All lanes RMDN1 pAb at 0.6ug/ml
Lane 1: K562 whole cell lysate
Lane 2: HepG-2 whole cell lysate
Lane 3: MCF7 whole cell lysate
Secondary: Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 36 kDa
Observed band size: 36 kDa



Immunohistochemistry of paraffin-embedded human skeletal muscle tissue using RMDN1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer using RMDN1 Polyclonal Antibody at dilution of 1:100