
MAMDC2 Polyclonal Antibody

(Catalog # A59746)

Background

endoplasmic reticulum

Description

MAMDC2 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

Q7Z304

Purification

>95%, Protein G purified

Immunogen

Recombinant Human MAM domain-containing protein 2 protein (455-686AA)

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

MAM domain-containing protein 2, MAMDC2, MAM domain-containing proteoglycan, Mamcan

Application

ELISA; Recommended dilution: ELISA 1:500-1:2000
