
MFGE8 Polyclonal Antibody, Biotin Conjugated

(Catalog # A53059)

Background

Plays an important role in the maintenance of intestinal epithelial homeostasis and the promotion of mucosal healing. Promotes VEGF-dependent neovascularization By similarity. Contributes to phagocytic removal of apoptotic cells in many tissues. Specific ligand for the alpha-v/beta-3 and alpha-v/beta-5 receptors. Also binds to phosphatidylserine-enriched cell surfaces in a receptor-independent manner. Zona pellucida-binding protein which may play a role in gamete interaction. Binds specifically to rotavirus and inhibits its replication. Ref. 15 Medin is the main constituent of aortic medial amyloid.

Description

MFGE8 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q08431

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Lactadherin protein (25-334AA)

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

Breast epithelial antigen BA46 HMFG MFGM Milk fat globule-EGF factor 8 Short name=MFG-E8 SED1 Cleaved into the following 2 chains: Lactadherin short form Medin

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
