
TL1A Monoclonal Antibody [L4G9]

(Catalog No. A-0608)

Background

TL1A is a cytokine that belongs to the tumor necrosis factor (TNF) ligand family. This protein is not expressed in B or T cells, but is liberally expressed in endothelial cells. The expression of this protein is brought on by TNF and IL-1 alpha. TL1A is a receptor for TNFRSF25 and TNFRSF6B. It can initiate NF-kappaB and MAP kinases, and acts as an autocrine factor to induce apoptosis in endothelial cells. TL1A also inhibits endothelial cell proliferation, and may function as an angiogenesis inhibitor. Two transcript variants encoding different isoforms have been found for this gene. Diseases associated with TL1A include primary biliary cirrhosis and diverticulitis.

Description

Mouse monoclonal antibody to TL1A, Clone L4G9

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant human TL1A extracellular domain.

Uniprot ID

O95150

Specificity

Human

Isotype

IgG1

Purification

Protein G- affinity purified

Storage

Store at 4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

Application

ELISA, FC (1:1000)

Ordering Information

Products

TL1A Monoclonal Antibody [L4G9]

Size

100 µg

Cat. No.

A-0608-100

This product is for research purposes only. Not intended for use in diagnostic procedures.