
TREML2 Antibody

(Catalog #A61486)

Background

Cell surface receptor that may play a role in the innate and adaptive immune response. Acts as a counter-receptor for CD276 and interaction with CD276 on T-cells enhances T-cell activation.

Description

TREML2 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

Q5T2D2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Trem-like transcript 2 protein (19-268AA)

Storage

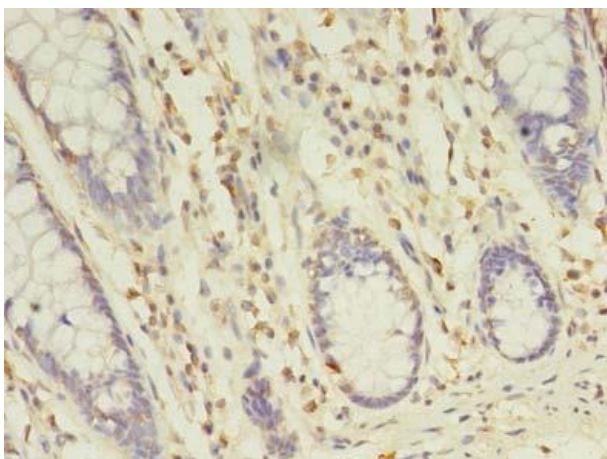
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

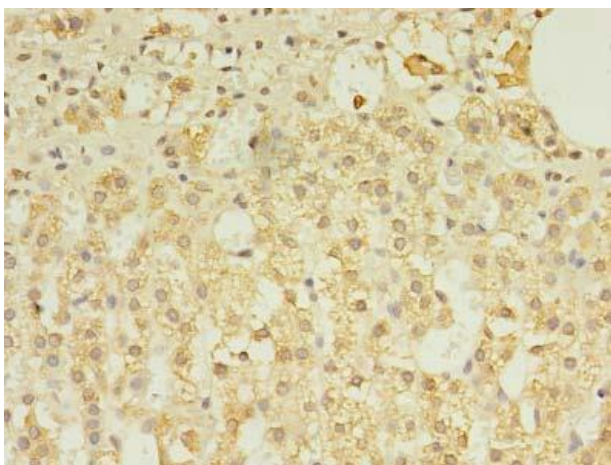
Trem-like transcript 2 protein, TLT-2, TREML2, C6orf76, TLT2, UNQ6268/PRO20473, Triggering receptor expressed on myeloid cells-like protein 2

Application

ELISA, IHC; Recommended dilution: IHC:1:20 - 1:200



Immunohistochemistry of paraffin-embedded human colon cancer using TREML2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using TREML2 Polyclonal Antibody at dilution of 1:100