

TRBC1 Polyclonal Antibody

(Catalog # A56723)

TRBC1 (T Cell Receptor Beta Constant 1) is a protein coding gene. Among its related pathways are Immunoregulatory interactions between a Lymphoid and a non-Lymphoid cell and Translocation of ZAP-70 to Immunological synapse. An important paralog of this gene is TRBC2.

Description

TRBC1 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

P01850

Purification

Protein G purified

Immunogen

Recombinant Human T cell receptor beta constant 1 protein (1-176AA)

Storage

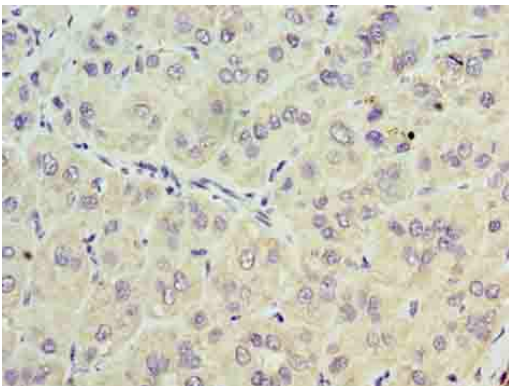
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

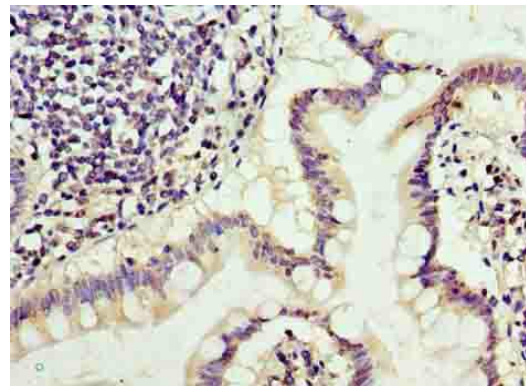
TRBC1

Application

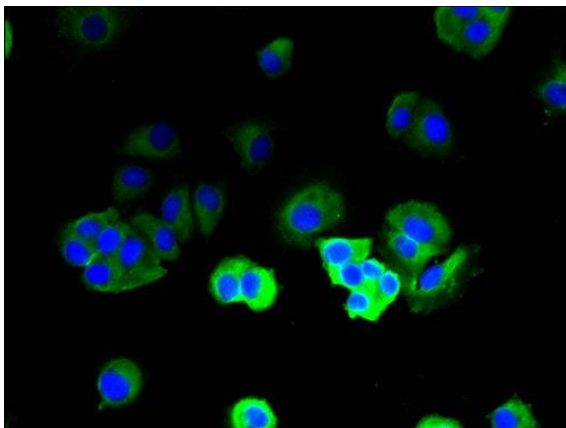
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human liver cancer at dilution of 1:100



Immunohistochemistry of paraffin-embedded human small intestine tissue at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)