

TIGIT Polyclonal Antibody

(Catalog # A64998)

Background

Binds with high affinity to the poliovirus receptor (PVR) which causes increased secretion of IL10 and decreased secretion of IL12B and suppresses T-cell activation by promoting the generation of mature immunoregulatory dendritic cells.

Description

TIGIT 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

Q495A1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human T-cell immunoreceptor with Ig and ITIM domains protein (22-141AA)

Storage

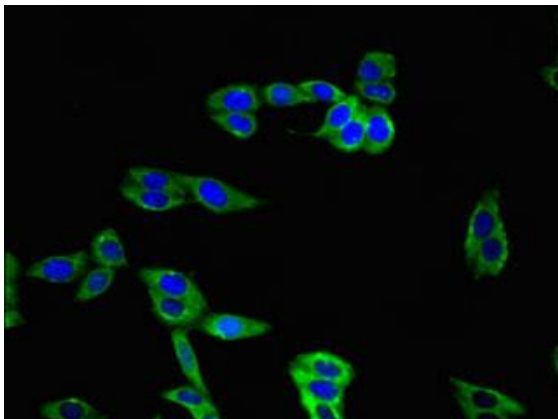
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

T-cell immunoreceptor with Ig and ITIM domains, V-set and immunoglobulin domain-containing protein 9, V-set and transmembrane domain-containing protein 3, TIGIT, VSIG9, VSTM3

Application

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



Immunofluorescent analysis of HepG2 cells using TIGIT Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)