

PVRIG Polyclonal Antibody

(Catalog # A65226)

Background

PVRIG (PVR Related Immunoglobulin Domain Containing) is a cell surface receptor for NECTIN2. It may act as a coinhibitory receptor that suppresses T-cell receptor-mediated signals. Following interaction with NECTIN2, this gene inhibits T-cell proliferation. It competes with CD226 for NECTIN2-binding.

Description

PVRIG Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q6DKI7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane protein PVRIG protein (193-326AA)

Storage

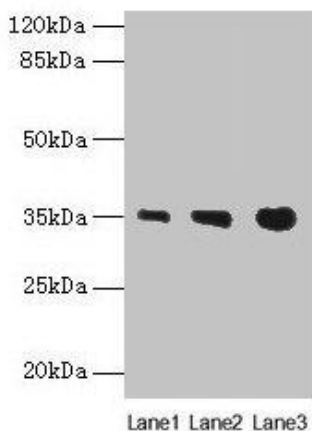
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transmembrane protein PVRIG, Poliovirus receptor-related immunoglobulin domain-containing protein, PVRIG, C7orf15

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



Western blot

All lanes: PVRIG antibody at 4μg/ml

Lane 1: Mouse liver tissue

Lane 2: Mouse kidney tissue

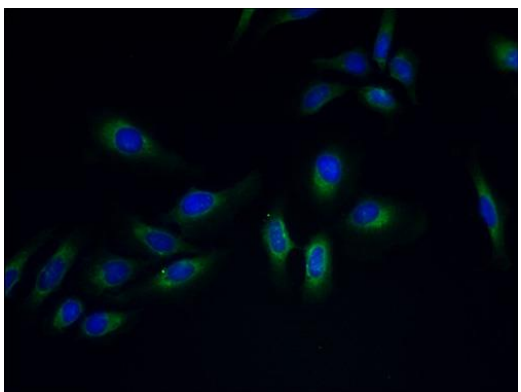
Lane 3: Mouse heart tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa



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