

PILRA Polyclonal Antibody

(Catalog # A68106)

Background

Paired receptors consist of highly related activating and inhibitory receptors and are widely involved in the regulation of the immune system. PILRA is thought to act as a cellular signaling inhibitory receptor by recruiting cytoplasmic phosphatases like PTPN6/SHP-1 and PTPN11/SHP-2 via their SH2 domains that block signal transduction through dephosphorylation of signaling molecules. Receptor for PIANP.

Description

PILRA 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

Q9UKJ1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Paired immunoglobulin-like type 2 receptor alpha protein (22-226AA)

Storage

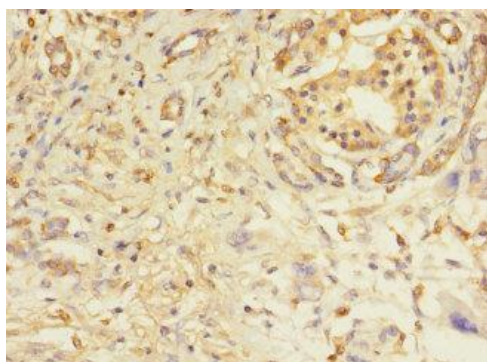
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Paired immunoglobulin-like type 2 receptor alpha, Cell surface receptor FDF03, Inhibitory receptor PILR-alpha, PILRA

Application

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



Immunohistochemistry of paraffin-embedded human pancreatic tissue PILRA Polyclonal Antibody at dilution 1:100