

LAIR2 Polyclonal Antibody

(Catalog #A62862)

Background

extracellular region, regulation of immune response

Description

LAIR2 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

Q6ISS4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Leukocyte-associated immunoglobulin-like receptor 2 protein (22-152AA)

Storage

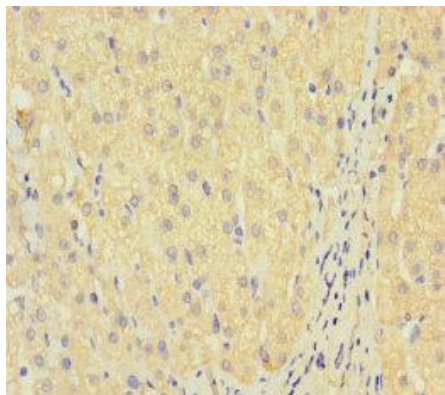
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Leukocyte-associated immunoglobulin-like receptor 2, LAIR-2, CD306, LAIR2, CD306

Application

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



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