
KIR2DS1 Polyclonal Antibody

(Catalog #A57497)

Background

Receptor on natural killer (NK) cells for HLA-C alleles. Does not inhibit the activity of NK cells.

Description

KIR2DS1 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

Q14954

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Killer cell immunoglobulin-like receptor 2DS1 protein (22-245AA)

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

CD158 antigen-like family member H, MHC class I NK cell receptor Eb6 Actl, CD158h, KIR2DS1

Application

ELISA; Recommended Dilution: ELISA 1:2000 – 1:5000
