

HLA-G Polyclonal Antibody

(Catalog # A53328)

Background

HLA class I histocompatibility antigen, alpha chain G (HLA-G) is involved in the presentation of foreign antigens to the immune system. It plays a role in maternal tolerance of the fetus by mediating protection from the deleterious effects of natural killer cells, cytotoxic T-lymphocytes, macrophages and mononuclear cells. Diseases associated with HLA-G include asthma susceptibility, Hla-G-related and severe pre-eclampsia. Among its related pathways are Interferon gamma signaling and HTLV-I infection.

Description

HLA-G 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

P17693

Purification

>95%, Protein G purified

Immunogen

Recombinant Human HLA class I histocompatibility antigen, alpha chain G protein (25-337AA)

Storage

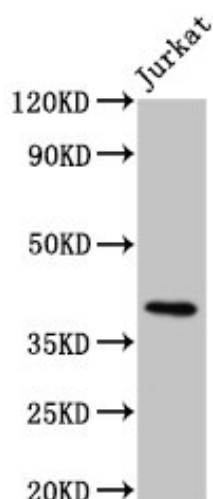
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

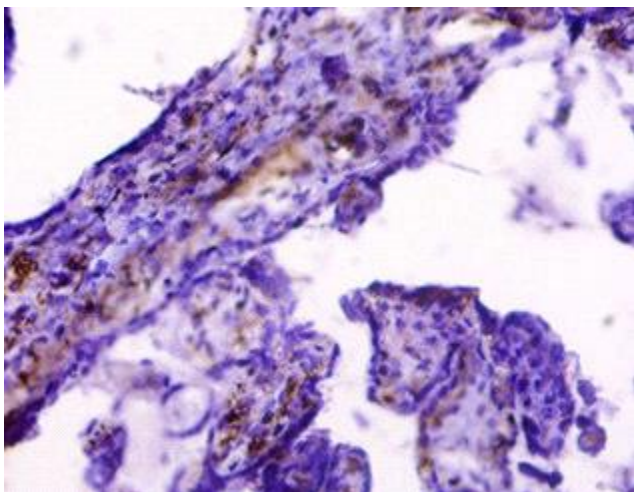
HLA G antigen, MHC class I antigen G, HLA-G, HLA-6.0, HLAG

Application

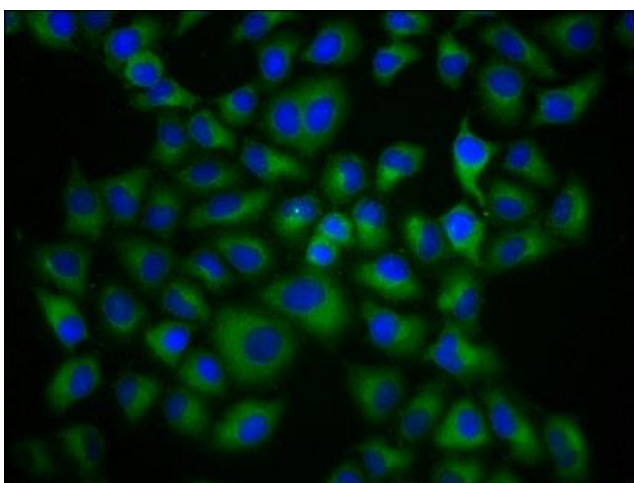
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-2000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot
Positive WB detected in: Jurkat whole cell lysate
All lanes: HLA-G polyclonal antibody at 8µg/ml
Secondary
Goat polyclonal to rabbit IgG at 1/50000 dilution
Predicted band size: 39 kDa
Observed band size: 39 kDa



Immunohistochemistry of paraffin-embedded human placenta tissue using HLA-G Polyclonal Antibody at dilution of 1:50.



Immunofluorescence staining of HeLa cells with HLA-G Polyclonal Antibody at 1:133, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).