
FCGRT Polyclonal Antibody

(Catalog # A71258)

Background

This gene encodes a receptor that binds the Fc region of monomeric immunoglobulin G. The encoded protein transfers immunoglobulin G antibodies from mother to fetus across the placenta. This protein also binds immunoglobulin G to protect the antibody from degradation. Alternative splicing results in multiple transcript variants.

Description

FCGRT Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P55899

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 25-300 of human FCGRT (NP_004098.1).

Storage

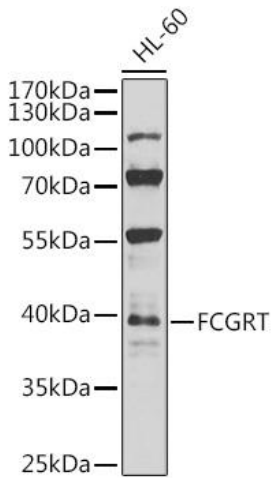
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

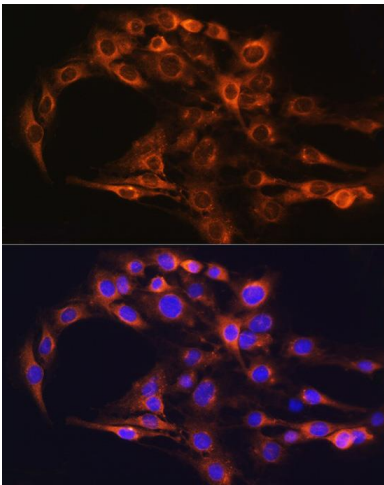
FCGRT; FCRN; alpha-chain; IgG receptor FcRn large subunit p51

Application

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



Western blot analysis of extracts of HL-60 cells, using FCGRT Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of C6 cells using FCGRT Polyclonal Rabbit pAb (A8544) at dilution of 1:100 (40x lens).
Blue: DAPI for nuclear staining.