

## FCGRT Polyclonal Antibody

(Catalog # A53808)

### Background

Binds to the Fc region of monomeric immunoglobulins gamma. Mediates the selective uptake of IgG from milk and helps newborn animals to acquire passive immunity. IgG in the milk is bound at the apical surface of the intestinal epithelium. The resultant FcRn-IgG complexes are transcytosed across the intestinal epithelium and IgG is released from FcRn into blood or tissue fluids (By similarity). Possible role in transfer of immunoglobulin G from mother to fetus.

### Description

FCGRT Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: Liquid in PBS containing 50% glycerol, 0.5% BSA\* and 0.02% sodium azide

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

P55899

### Purification

Antigen affinity purification

### Immunogen

Synthesized peptide derived from human FCGRN

### Storage

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

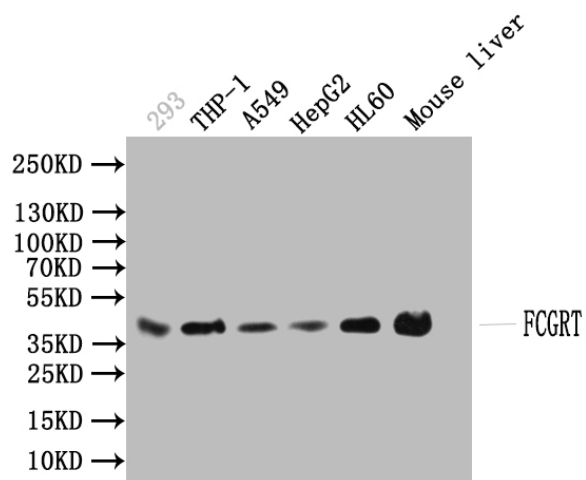
### Alternative Names

IgG Fc fragment receptor transporter alpha chain FCGRT FCRN

### Application

ELISA, WB; Recommended dilution: WB:1:1000-1:3000

*\*The BSA contained in this antibody is exempt from the requirement of certification by an authoritative body. This exemption is due to the comprehensive purification process, which ensures the complete absence of viable microorganisms, the BSA's sourcing from aseptically collected serum in the USA, and its subsequent sterile filtration and lyophilization. Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*



#### Western Blot

Positive WB detected in: 293 whole cell lysate, THP-1 whole cell lysate, A549 whole cell lysate, HepG2 whole cell lysate, HL60 whole cell lysate, Mouse Liver tissue lysate

All lanes: FCGRT Polyclonal Antibody at 1:1000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 40kDa

Observed band size: 40kDa