

---

## KLRG1 Polyclonal Antibody, Biotin Conjugated

(Catalog # A55658)

---

### Background

Plays an inhibitory role on natural killer (NK) cells and T-cell functions upon binding to their non-MHC ligands. May mediate missing self-recognition by binding to a highly conserved site on classical cadherins, enabling it to monitor expression of E-cadherin/CDH1, N-cadherin/CDH2 and R-cadherin/CDH4 on target cells.

### Description

KLRG1 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q96E93

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Killer cell lectin-like receptor subfamily G member 1 protein (60-195AA)

### 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

C-type lectin domain family 15 member A, ITIM-containing receptor MAFA-L, MAFA-like receptor, Mast cell function-associated antigen, KLRG1, CLEC15A, MAFA, MAFAL

### Application

ELISA; Recommended dilution: ELISA 1:500-1:1000

---