
LY6G6D Recombinant Monoclonal Antibody [8I8]

(Catalog #A74075)

Background

LY6G6D belongs to a cluster of leukocyte antigen-6 (LY6) genes located in the major histocompatibility complex (MHC) class III region on chromosome 6. Members of the LY6 superfamily typically contain 70 to 80 amino acids, including 8 to 10 cysteines. Most LY6 proteins are attached to the cell surface by a glycosylphosphatidylinositol (GPI) anchor that is directly involved in signal transduction.

Description

LY6G6D Recombinant Monoclonal Antibody [8I8]. Unconjugated. Raised in: HEK293F Cell.

Formulation

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

Specificity

Human, Macaca fascicularis

Isotype

hIgG1

Uniprot ID

O95868

Purification

Affinity-chromatography

Immunogen

Recombinant Human LY6G6D protein

Storage

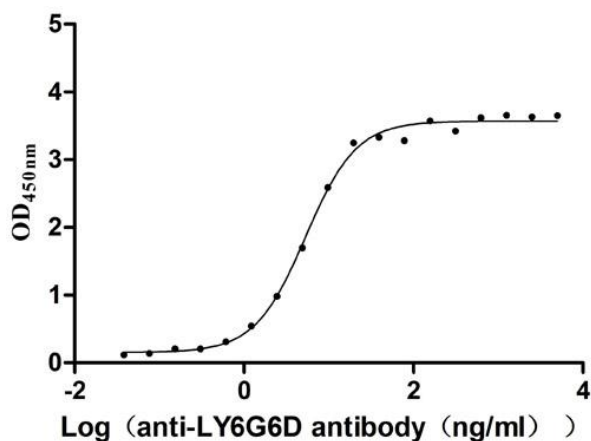
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

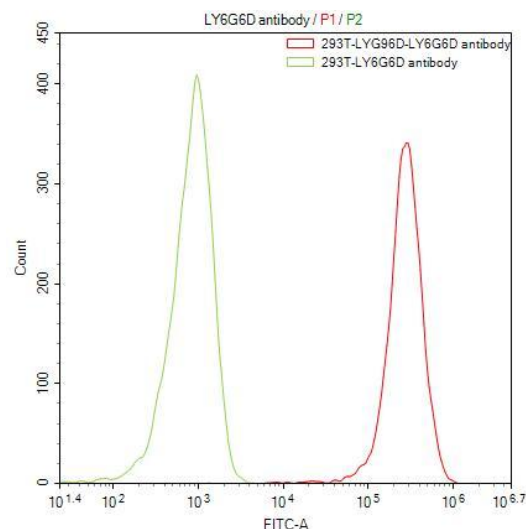
Lymphocyte antigen 6 complex locus protein G6d; Protein Ly6-D; Megakaryocyte-enhanced gene transcript 1 protein; LY6G6D; C6orf23; G6D; MEGT1; NG25

Application

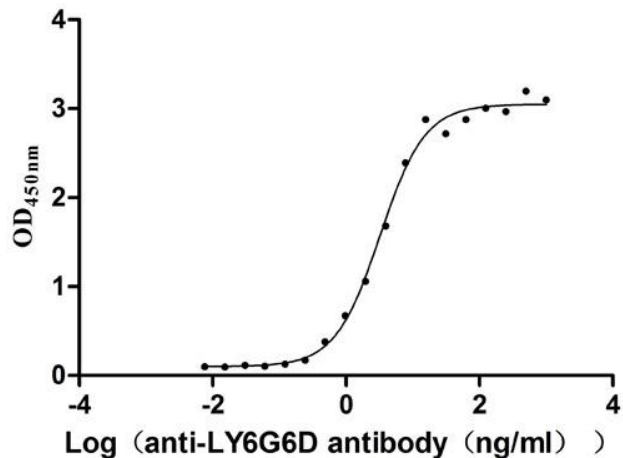
ELISA, FC; Recommended dilution: ELISA:1:500-1:5000, FC:1:100 – 1:1000



The Binding Activity of LY6G6D with anti-LY6G6D antibody Activity: Measured by its binding ability in a functional ELISA. Immobilized Macaca fascicularis LY6G6D at 2 ug/mL can bind Anti-LY6G6D recombinant antibody, the EC₅₀ is 3.963-7.154 ng/mL.



Untransfected HEK293T cells (green line) and transfected Human LY6G6D HEK293T stable cells (red line) were stained with anti-LY6G6D antibody (LY6G6D rMAb) (2μg/1*10⁶cells), washed and then followed by FITC-conjugated anti-Human IgG Fc antibody and analyzed with flow cytometry.



The Binding Activity of LY6G6D with anti-LY6G6D antibody Activity: Measured by its binding ability in a functional ELISA. Immobilized Human LY6G6D at 2 ug/mL can bind Anti-LY6G6D recombinant antibody, the EC₅₀ is 2.816-3.741 ng/mL.