
FCER1G Polyclonal Antibody, Biotin Conjugated

(Catalog #A54648)

Background

Associates with a variety of FcR alpha chains to form a functional signaling complex. Regulates several aspects of the immune response. The gamma subunit has a critical role in allowing the IgE Fc receptor to reach the cell surface. Also involved in collagen-mediated platelet activation and in neutrophil activation mediated by integrin.

Description

FCER1G 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

P30273

Purification

Protein G purified

Immunogen

Recombinant Human High affinity immunoglobulin epsilon receptor subunit gamma protein (19-86AA)

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

Fc receptor gamma-chain Fc-epsilon RI-gamma IgE Fc receptor subunit gamma FCER1G

Application

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