
NKG7 Polyclonal Antibody, Biotin Conjugated

(Catalog #A60027)

Background

integral component of plasma membrane

Description

NKG7 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

Q16617

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein NKG7 protein (30-60AA)

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

Protein NKG7, NKG7, GIG1, G-CSF-induced gene 1 protein, GIG-1 protein, Granule membrane protein of 17 kDa, GMP-17, Natural killer cell protein 7, p15-TIA-1

Application

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