
Cd63 Polyclonal Antibody, HRP Conjugated

(Catalog # A53108)

Background

May play some role in signal transduction pathways.

Description

Cd63 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P41731

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse CD63 antigen protein (103-203AA)

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

CD_antigen=CD63, Cd63

Application

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