

H2-D1 Polyclonal Antibody

(Catalog #A57413)

Background

Involved in the presentation of foreign antigens to the immune system.

Description

H2-D1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P01900

Purification

Protein G purified

Immunogen

Recombinant Mouse H-2 class I histocompatibility antigen, D-D alpha chain protein (25-311AA)

Storage

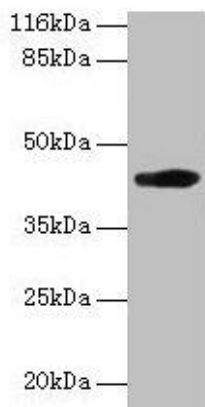
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

H-2D(D), H2-D1

Application

ELISA, WB; Recommended dilution: WB:1:500-1:5000



Western blot

All lanes: H2-D1 antibody at 2ug/ml + A549 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 42 kDa

Observed band size: 42 kDa