
H2-K1 Polyclonal Antibody

(Catalog #A53742)

Description

H2-K1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P01902

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse H-2 class I histocompatibility antigen, K-D alpha chain protein (22-305AA)

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

H2-K1 , H2-K

Application

ELISA; Recommended Dilution: ELISA 1:2000 – 1:5000
