
HLA-B Polyclonal Antibody, FITC Conjugated

(Catalog # A52235)

Background

Involved in the presentation of foreign antigens to the immune system

Description

HLA-B Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q96DW9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human HLA-B protein (25-308AA)

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

HLA-B

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
