
HLA-E Polyclonal Antibody, FITC Conjugated

(Catalog # A50750)

Description

HLA-E 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

P13747

Purification

Affinity purified

Immunogen

Recombinant Human HLA class I histocompatibility antigen, alpha chain E protein (33-261AA)

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 class I histocompatibility antigen, alpha chain E, MHC class I antigen E, HLA-E, HLA-6.2, HLAE

Application

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