

ADA Polyclonal Antibody

(Catalog # A70438)

Background

This gene encodes an enzyme that catalyzes the hydrolysis of adenosine to inosine. Various mutations have been described for this gene and have been linked to human diseases. Deficiency in this enzyme causes a form of severe combined immunodeficiency disease (SCID), in which there is dysfunction of both B and T lymphocytes with impaired cellular immunity and decreased production of immunoglobulins, whereas elevated levels of this enzyme have been associated with congenital hemolytic anemia.

Description

ADA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P00813

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-363 of human ADA (NP_000013.2).

Storage

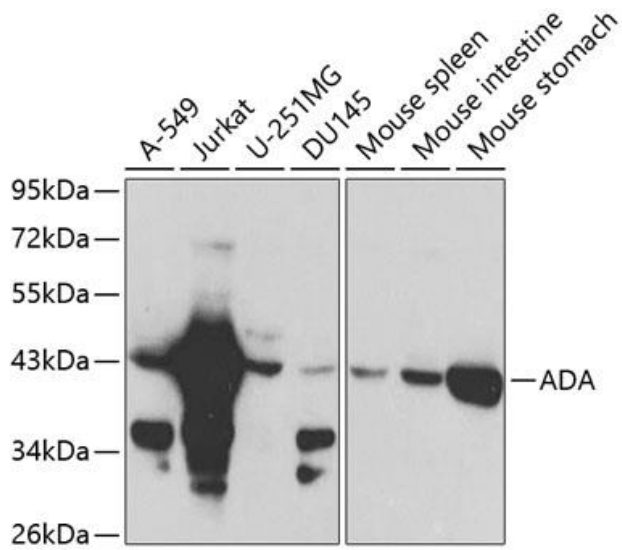
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

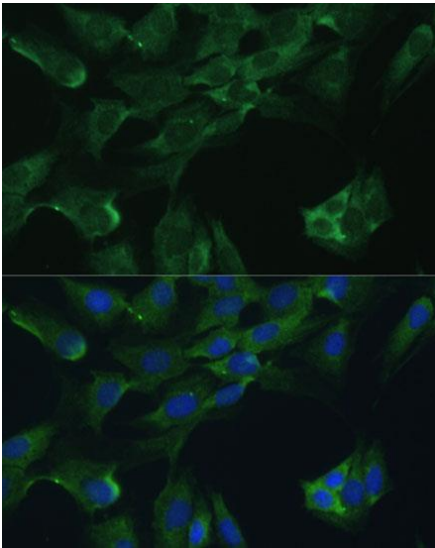
ADA; adenosine deaminase

Application

WB, IF; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200



Western blot analysis of extracts of various cell lines, using ADA Polyclonal Antibody at 1:1000 dilution.
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of C6 cells using ADA Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.