
IL2 Polyclonal Antibody, FITC Conjugated

(Catalog # A59476)

Background

Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

Description

IL2 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

P60568

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Interleukin-2 protein (21-153AA)

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

Interleukin-2, IL-2, IL2, T-cell growth factor, TCGF, INN: Aldesleukin

Application

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