
IL4 Polyclonal Antibody

(Catalog # A70790)

Background

The protein encoded by this gene is a pleiotropic cytokine produced by activated T cells. This cytokine is a ligand for interleukin 4 receptor. The interleukin 4 receptor also binds to IL13, which may contribute to many overlapping functions of this cytokine and IL13. STAT6, a signal transducer and activator of transcription, has been shown to play a central role in mediating the immune regulatory signal of this cytokine. This gene, IL3, IL5, IL13, and CSF2 form a cytokine gene cluster on chromosome 5q, with this gene particularly close to IL13. This gene, IL13 and IL5 are found to be regulated coordinately by several long-range regulatory elements in an over 120 kilobase range on the chromosome. Two alternatively spliced transcript variants of this gene encoding distinct isoforms have been reported.

Description

IL4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P05112

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 25-153 of human IL4 (NP_000580.1).

Storage

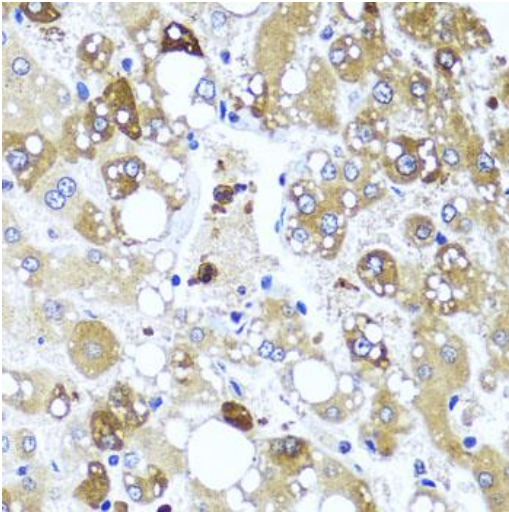
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

IL4; BCGF-1; BCGF1; BSF-1; BSF1; IL-4; interleukin-4

Application

WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100



Immunohistochemistry of paraffin-embedded human liver damage using IL4 Polyclonal Antibody at dilution of 1:100 (40x lens).