

Human IL-4 (recombinant)



Alias: none **Catalog #:** 6486

Size: 5 ug **Research Use Only**

Molecular Weight: 15.0 kDa

Source: Yeast. Recombinant Human IL-4 was produced in yeast and, therefore, does not have endotoxin. It is naturally folded and post-translationally modified.

Formulation: Lyophilized without carrier protein.

Purity: >95% as visualized by SDS-PAGE analysis.

Purification: Ion-exchange chromatography.

Bioactivity: In testing.

Entrez Gene ID: 3565

Number of Amino Acids: 129

Amino Acid Sequence: HKCDITLQEI IKTLNSLTEQ KTLCTELTVT DIFAASKNTT EKETFCRAAT VLRQFYSHHE
KDTRCLGATA QQFHRHKQLI RFLKRLDRNL WGLAGLNSCP VKEANQSTLE NFLERLKTIM
REKYSKCSS (129)

Country of Origin: USA

Reconstitution: Reconstitute with sterile phosphate-buffered saline containing at least 0.1% carrier protein.

Stability and Storage: Stable for up to twelve months from date of receipt at -20°C. Stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C. Avoid repeated freeze/thaw cycles.

Applications: The Human IL-4 protein can be used in cell culture, as an IL-4 ELISA Standard, and as a Western Blot Control.

Background: Interleukin-4 (IL-4) induces differentiation of naive helper T cells (Th0 cells) to Th2 cells. Upon activation by IL-4, Th2 cells subsequently produce additional IL-4. IL-4 has many biological roles, including the stimulation of activated B-cell and T-cell proliferation, and the differentiation of CD4+ T-cells into Th2 cells. It is a key regulator in humoral and adaptive immunity. IL-4 induces B-cell class switching to IgE, and up-regulates MHC class II production. Overproduction of IL-4 is associated with allergies.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Human IL-4 updated on: 1/29/2014