IL-4 (mouse), (recombinant)

Interleukin 4 is a cytokine that has various biological functions including stimulation of activated B-cells, stimulation of T-cell proliferation, and the differentiation of B cells into plasma cells. It is a key regulator in humoral and adaptive immunity. IL-4 decreases the production of Th1 cells, macrophages, IFN- γ , and dendritic cell IL-12. Overexpression of IL-4 is associated with allergies.

Citations: 2

View Online »

Ordering Information

Order Online »

ENZ-PRT133-0020

20µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-4

Formulation Lyophilized.

MW ~14 kDa

Purity ≥98% (SDS-PAGE gel & HPLC)

Produced in E.coli. Source

UniProt ID P07750

Last modified: May 29, 2024



info-

eu@enzolifesciences.com