

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	20 μ g
-----------------	------------

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)

Source Produced in *E.coli*.

UniProt ID P07750

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)