

IL-10 (human), (recombinant)

IL-10, also known as human cytokine synthesis inhibitory factor (CSIF), is an anti-inflammatory cytokine. IL-10 inhibits the expression of proinflammatory cytokines such as IL-1 and TNF- α . Like IL-4, IL-10 enhances humoral immune responses and attenuates cell-mediated immune reactions.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT134-0010	10 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-10, Cytokine synthesis inhibitory factor, CSIF

Formulation Lyophilized from 5mM sodium phosphate, pH 7.2.

MW ~18.6 kDa

Purity ≥98% (SDS-PAGE gel & HPLC)

Source Produced in *E.coli*.

UniProt ID P22301

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)