

IL-11 (mouse), (recombinant)

Interleukin 11 (IL-11) is a member of the gp130 family of cytokines. IL-11 functions to promote hematopoietic stem cell proliferation and megakaryocyte differentiation. IL-11 binds the IL-11R α receptor to activate JAK downstream signaling. In non-hematopoietic cell populations, IL-11 stimulates acute-phase proteins, modulates the development of immunoglobulin-producing B cells, and regulates bone turnover.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

 RUO - Research Use Only

Product Details

Alternative Name	Interleukin-11, AGIF, Adipogenesis Inhibitory Factor, Oprelvekin
Biological Activity	T11 cell proliferation; ED50 \leq 2.5 ng/mL (\geq 4.0 x 10 ⁵ units/mg).
Endotoxin Content	\leq 1 EU/ μ g protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 0.1% Trifluoroacetic Acid (TFA).
MW	~19.3 kDa
Purity	\geq 95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 179 amino acids.
UniProt ID	P47873

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)