

IL-1 β (human), (recombinant) (cell culture grade)

Potent immunomodulator

Potent immunomodulator. Mediates a wide range of immune and inflammatory responses including the activation of B and T cells.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-520-0010	10 μ g
--------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	IL-1F2, Interleukin-1 β
Biological Activity	Exerts <i>in vitro</i> biological activity in the range of 0.1-10ng/ml.
CAS	97599-21-8
Formulation	Lyophilized. Salt-free.
MI	14: 4993
MW	~17kDa.
Purity	≥97% (SDS-PAGE)
Reconstitution	Reconstitute with water, PBS or aqueous buffer to a concentration of 100 μ g/ml. For further dilution carrier protein (~1%) should be added to avoid loss of bioactivity.
Source	Produced in <i>E. coli</i> .
Specific Activity	≥1×10 ⁷ U/mg protein. Determined by proliferation assay using human TF-1 (T. Kitamura, et al.; Int. Immunol. 3 , 571 (1991)). The ED ₅₀ for this effect is typically 0.2-2ng/ml. Results were compared with the international IL-1 reference standard preparation (WHO 86/680, NIBSC).

UniProt ID P01584



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

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

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

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

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

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

Last modified: May 29, 2024