

# IL-1 $\beta$ (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 $\mu$ 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 $\beta$
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	<b>14:</b> 4993
MW	~17kDa.
Purity	≥97% (SDS-PAGE)
Reconstitution	Reconstitute with water, PBS or aqueous buffer to a concentration of 100 $\mu$ 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 <sup>7</sup> U/mg protein. Determined by proliferation assay using human TF-1 (T. Kitamura, et al.; Int. Immunol. <b>3</b> , 571 (1991)). The ED <sub>50</sub> 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](mailto: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](mailto:eu@enzolifesciences.com)

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

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

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

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

Last modified: May 29, 2024