

# IL-1 $\alpha$ (mouse), (recombinant)

Interleukin-1 $\alpha$  (IL-1 $\alpha$ ) is a protein of the interleukin-1 family. In general, Interleukin 1 is responsible for the production of inflammation, as well as the promotion of fever and sepsis. IL-1 $\alpha$  inhibitors are being developed to interrupt those processes and treat diseases.

**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-1 $\alpha$ , Hematopoietin-1, Lymphocyte-activating factor, LAF, Endogenous pyrogen, Leukocyte endogenous mediator, LEM, Mononuclear cell factor, MCF, IL-1F1
Formulation	Lyophilized from 1mM TRIS, pH 8.0.
MW	~17.9 kDa
Purity	≥98% (SDS & HPLC)
Purity Detail	Sterile filtered through a 0.2 micron filter.
Source	Produced in <i>E. coli</i> .
UniProt ID	P01582

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)