

IL-1 α (mouse), (recombinant)

Interleukin-1 α (IL-1 α) 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 α inhibitors are being developed to interrupt those processes and treat diseases.

Ordering Information

[Order Online »](#)

ENZ-PRT187-0010	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-1 α , 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)