# IL-1α (human),(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.

# **Ordering Information**

Order Online »

ENZ-PRT186-0010

10µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Handling** Centrifuge vial before opening. Do not vortex. For prolonged storage, dilute to working

aliquots in a 0.1% BSA solution, store at -20°C to -80°C. Avoid freeze/thaw cycles.

Short Term Storage -20°C

Long Term Storage -20°C

**Shipping** Ambient Temperature

### Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Interleukin-1α, Hematopoietin-1, Lymphocyte-activating

factor, LAF, Endogenous pyrogen, Leukocyte endogenous

mediator, LEM, Mononuclear cell factor, MCF, IL-1F1

Biological Activity Stimulates the proliferation of mouse D10S cells.

Formulation Lyophilized from 10mM Sodium Phosphate pH 8.0,

containing 50mM NaCl.

**MW** ~18.0 kDa

Purity ≥98% (SDS & HPLC)

Purity Detail Sterile filtered through a 0.2 micron filter.

**Reconstitution** Reconstitute in water to a concentration of 0.1 – 1.0

mg/mL.

Source Produced in E. coli.

**Specific Activity**  $1 \times 10^9$  units/mg. Determined by the ability of IL-1 $\alpha$  to

stimulate the proliferation of mouse D10S cells. The

expected ED<sub>50</sub> is typically ≤0.001 ng/mL.

UniProt ID P01583



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com