IL-1α (mouse), (recombinant)

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

Ordering Information

Order Online »

ENZ-PRT187-0010

10µg

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-1α, 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 *E. coli*.

UniProt ID P01582

Last modified: May 29, 2024



uk@enzolifesciences.com