IL-1β (mouse),(recombinant)

IL-1 β is produced by activated macrophages as a proprotein and cleaved by caspase-1. IL-1 β is an important mediator of the inflammatory response, and is involved in a variety of cellular activities including: cell proliferation, differentiation, and apoptosis. The induction of cyclooxygenase-2 (PTGS2/COX2) by IL-1 β in the CNS is found to contribute to inflammatory pain hypersensitivity.

Ordering Information

Order Online »

ENZ-PRT136-0010

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-1β, Catabolin

Formulation Lyophilized from 10 mM sodium citrate, pH 4.0.

MW ~17.5 kDa

Purity ≥98% (SDS-PAGE gel & HPLC)

Source Produced in E.coli.

UniProt ID P10749

Last modified: May 29, 2024



eu@enzolifesciences.com