

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

IL-1 $\beta$  is produced by activated macrophages as a proprotein and cleaved by caspase-1. IL-1 $\beta$  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 $\beta$  in the CNS is found to contribute to inflammatory pain hypersensitivity.

## Ordering Information

[Order Online »](#)

ENZ-PRT136-0010	10 $\mu$ 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 $\beta$ , Catabolin

Formulation Lyophilized from 10 mM sodium citrate, pH 4.0.

MW ~17.5 kDa

Purity  $\geq$ 98% (SDS-PAGE gel & HPLC)

Source Produced in *E.coli*.

UniProt ID P10749



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)