

IL-1 β (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.

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 β , 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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com