## IL-1β recombinant monoclonal antibody (IL1B/7049R)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

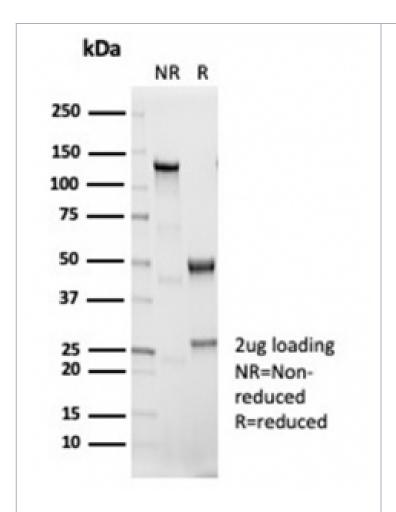
**Order Online** »

ENZ-ABS915-0100

100µg

Manuals, SDS & CofA

**View Online »** 



SDS-PAGE analysis of purified, BSA-free recombinant IL1 beta antibody (clone IL1B/7049R) as confirmation of integrity and purity.

## **Handling & Storage**

Handling Aliquot the recombinant IL1 beta antibody and store frozen at -20°C or colder. Avoid

freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Alternative Name** Interleukin-1β

**Application** ELISA, IHC (PS)

Clone IL1B/7049R

**Formulation** Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced), 0.05% sodium azide.

Host Rabbit

**Immunogen** Recombinant protein fragment of IL-1β.

Isotype IgG

**Purity Detail** Protein A/G affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (FFPE) (1-2µg/ml)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

**Species Reactivity** Human

**UniProt ID** P01584

This antibody is covered by our Worry-Free Guarantee **Worry-free Guarantee** 

Last modified: July 12, 2024



info-

eu@enzolifesciences.com