

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

This antibody is covered by our [Worry-Free Guarantee](#).

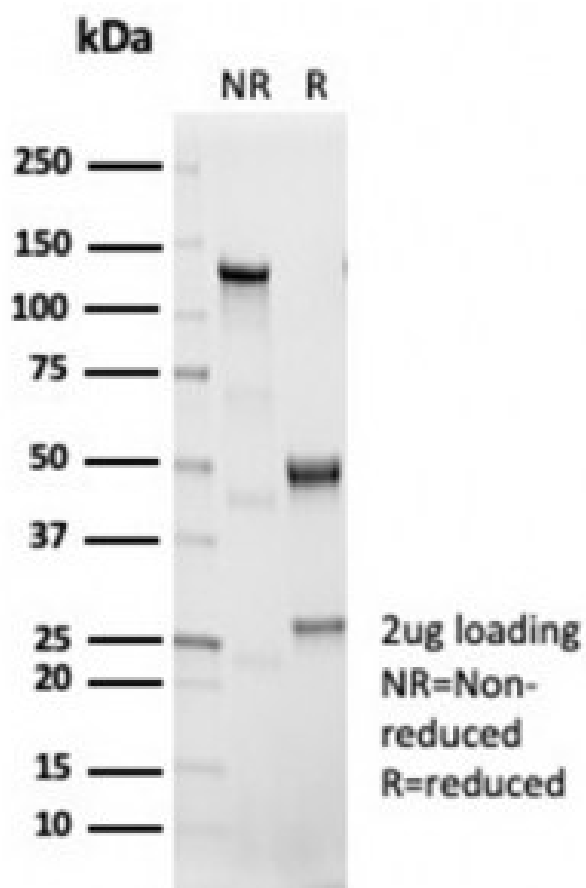
Ordering Information

[Order Online »](#)

ENZ-ABS915-0100	100 μ g
-----------------	-------------

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



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)