

# IL-21R (ED) polyclonal antibody

IL-21 is a cytokine identified in human and mouse that shows significant homology to IL-2, IL-4, and IL-15. The receptor for IL-21 (IL-21R) also termed NLR for novel interleukin receptor) is a member of the class I cytokine receptor family. IL-21R forms a complex with the common cytokine receptor gamma chain, gammac, and mediates IL-21 signaling via the JAK-Stat pathway. IL-21 is expressed in activated T-cells, as well as HL-60 and THP-1 cell lines. IL-21 plays a role in the proliferation and maturation of NK, B- and T-cell populations.

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

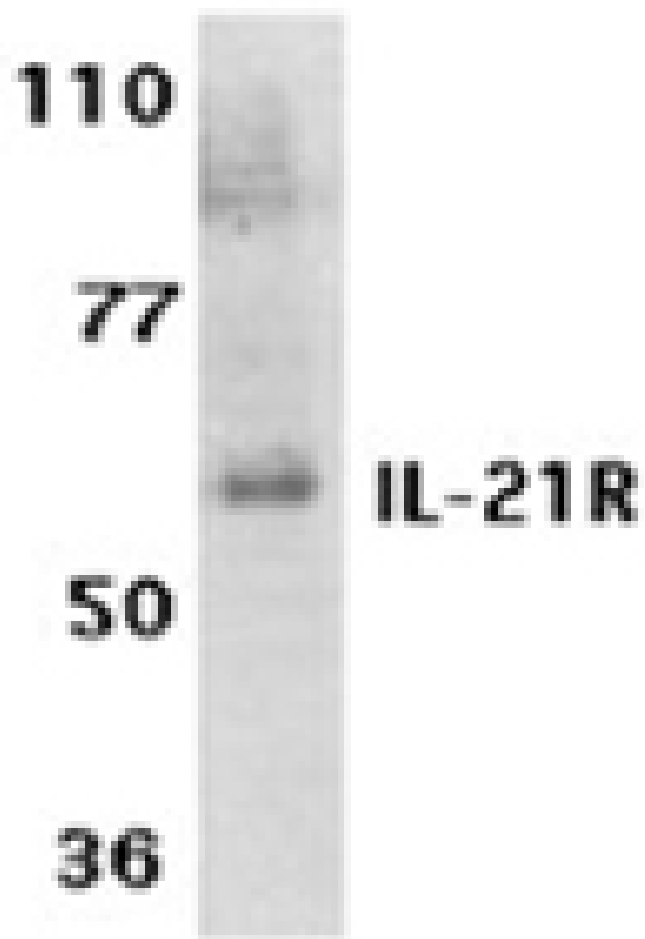
## Ordering Information

[Order Online »](#)

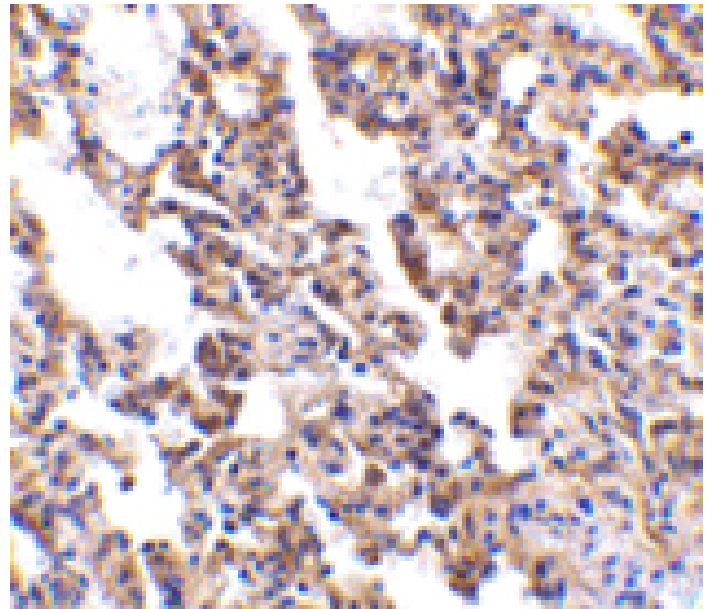
|                 |       |
|-----------------|-------|
| ADI-905-308-100 | 100µg |
|-----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of IL-21 receptor expression in human Raji cell lysate with IL-21 Receptor antibody at 1µg/ml.



Immunohistochemistry analysis of IL-21 receptor in rat lung with IL-21 receptor antibody at 10µg/ml.

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

|                                     |                                                                                                                                                                                                            |
|-------------------------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Alternative Name                    | Interleukin-21 receptor, IL-21 receptor, CD360                                                                                                                                                             |
| Application                         | IHC, WB                                                                                                                                                                                                    |
| Application Notes                   | Detects a band of ~60kDa by Western blot.                                                                                                                                                                  |
| Formulation                         | Liquid. In PBS containing 0.02% sodium azide.                                                                                                                                                              |
| Host                                | Rabbit                                                                                                                                                                                                     |
| Immunogen                           | Synthetic peptide corresponding to a portion of human IL-21R precursor.                                                                                                                                    |
| Purity Detail                       | Peptide affinity purified.                                                                                                                                                                                 |
| Recommendation Dilutions/Conditions | Immunohistochemistry (10µg/ml)Western Blot (0.5-1.0µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity                  | Human, Mouse, Rat                                                                                                                                                                                          |
| UniProt ID                          | Q9HBE5                                                                                                                                                                                                     |
| Worry-free Guarantee                | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>                                                                                                                                       |

Last modified: May 29, 2024



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-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@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)