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 NILR 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.

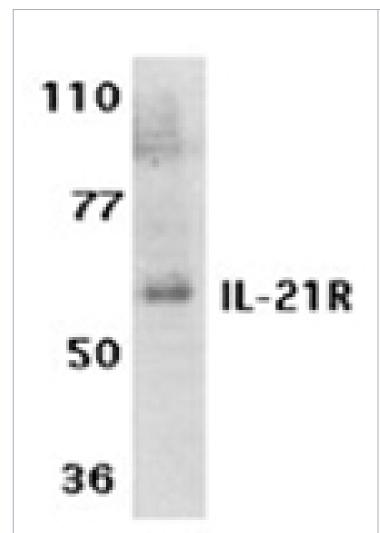
Ordering Information

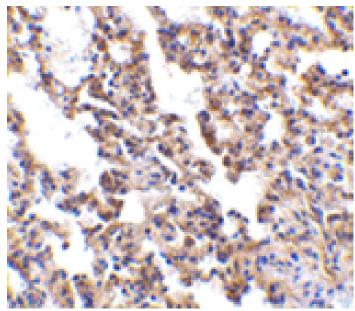
Order Online »

ADI-905-308-100

Manuals, SDS & CofA

View Online »





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

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

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 NotesDetects 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 Worry-Free Guarantee

•

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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