

IL-5RA polyclonal antibody

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

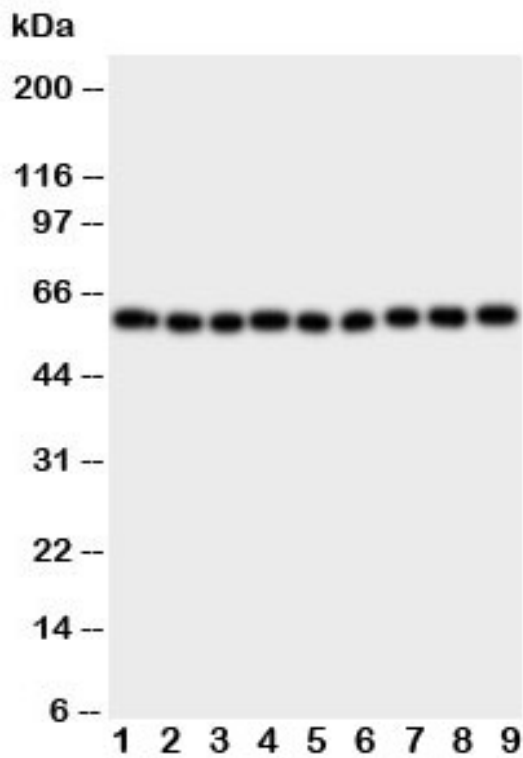
Ordering Information

[Order Online »](#)

ENZ-ABS900-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot testing of IL5RA antibody and Lane 1: A549; 2: SMMC-7721; 3: U87; 4: 293T; 5: PANC; 6: HeLa; 7: Jurkat; 8: Raji; 9: CEM cell lysate. The 420 amino acid protein is routinely visualized between 46~60KD, depending on level of glycosylation.

Handling & Storage

Use/Stability	After reconstitution, the IL5RA antibody can be stored for up to one month at +4°C. For long-term, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin-5 receptor subunit α , CD125
Application	WB
Formulation	Lyophilized from 1X PBS containing 2.5% BSA and 0.025% sodium azide/thimerosal.
Host	Rabbit
Immunogen	Amino acid sequence from the N-terminus of human IL-5RA (NPDQEQARNVNLEYQVK).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	Western blot (0.5-1 μ g/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	Q01344
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-
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)