

IL-4R polyclonal antibody

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

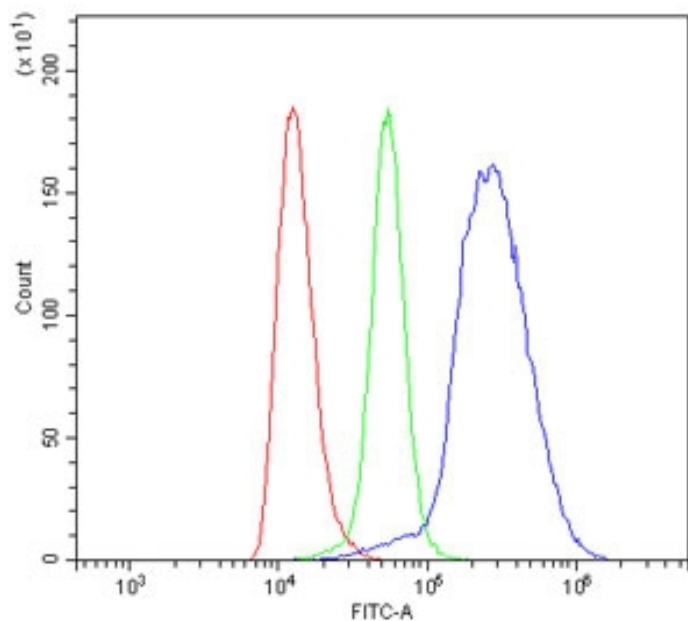
Ordering Information

[Order Online »](#)

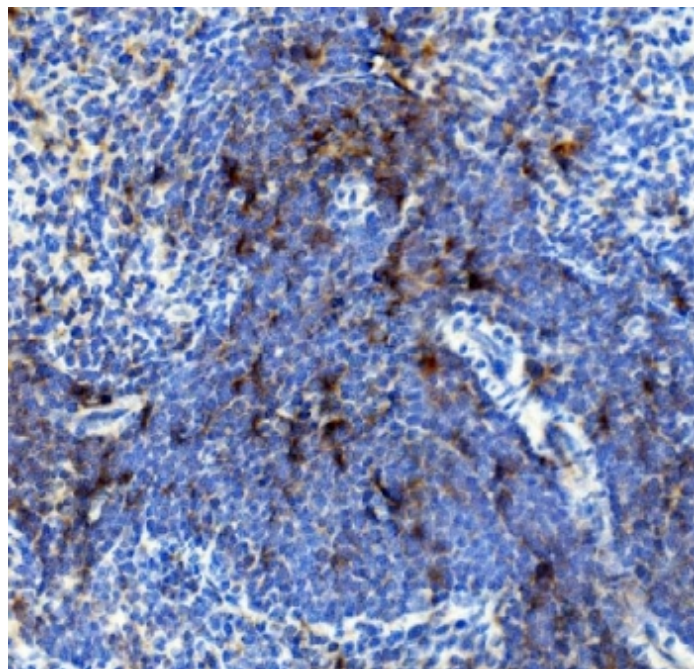
ENZ-ABS899-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry testing of mouse spleen cells with IL4 receptor antibody at $1\mu\text{g}/\text{million cells}$ (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= IL4 Receptor antibody.



IHC staining of FFPE rat spleen with IL4 receptor antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.

Handling & Storage

Use/Stability	After reconstitution, the IL4 receptor 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-4 receptor, CD124
Application	ELISA, Flow Cytometry, IHC (PS)
Formulation	Lyophilized from 1X PBS containing 2% Trehalose and 0.0125% sodium azide.
Host	Rabbit
Immunogen	Rat recombinant IL-4R (aa 26-719).
Isotype	IgG
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	ELISA (direct) (0.1-0.5µg/ml)Flow Cytometry (1-3µg/million cells)Immunohistochemistry (FFPE) (2-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse, Rat
UniProt ID	Q63257
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)