

IL-6 receptor recombinant monoclonal antibody (033)

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

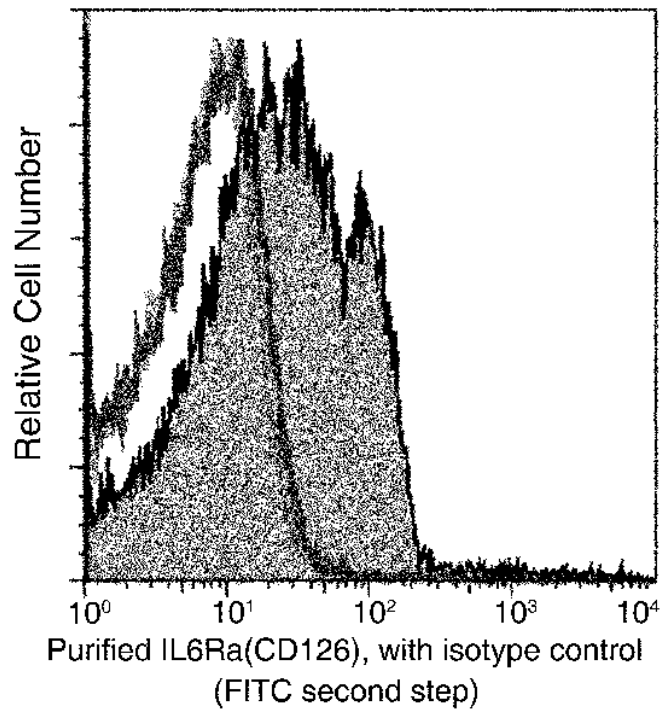
Ordering Information

[Order Online »](#)

ENZ-ABS1013-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse IL6Ra(CD126) expression on BABL/c splenocytes. Cells were stained with purified anti-Mouse IL6Ra(CD126), then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

Handling & Storage

Handling	Protected from light. Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CD126, IL-6R, Interleukin-6 receptor
Application	ELISA, Flow Cytometry
Clone	033
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant mouse IL-6R / CD126 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:25-1:100)ELISA (Capture) (1:250-1:2'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
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)