

# IL-17RB polyclonal antibody

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

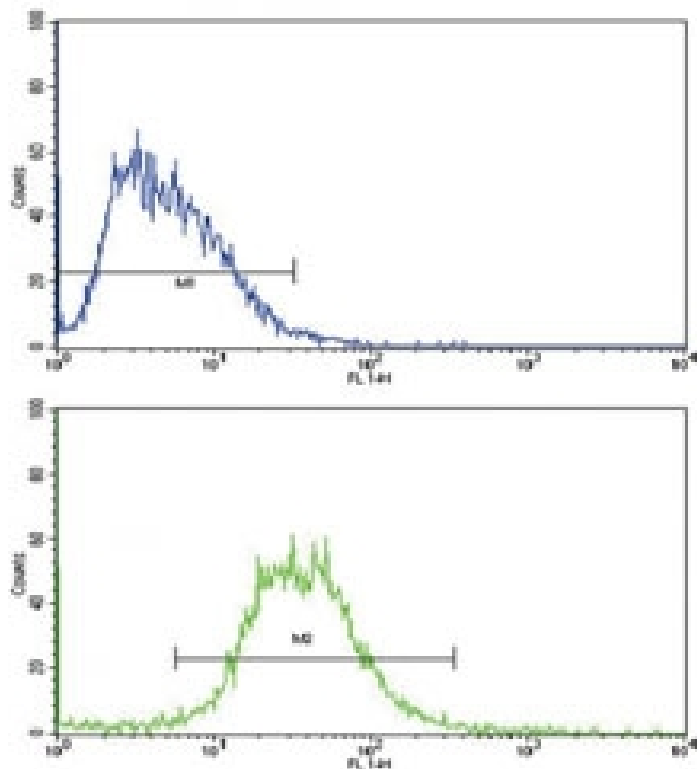
## Ordering Information

[Order Online »](#)

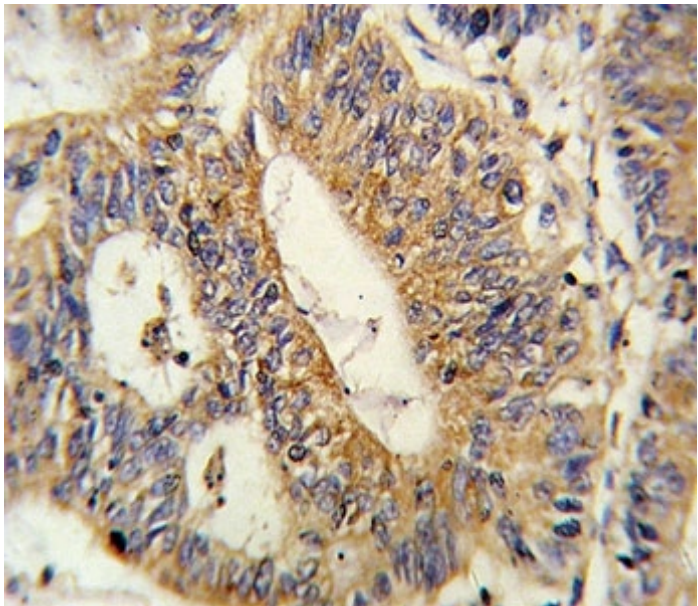
ENZ-ABS897-0400	0.4ml
-----------------	-------

## Manuals, SDS & CofA

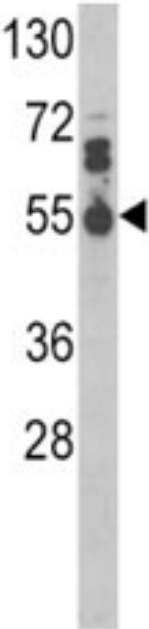
[View Online »](#)



Flow cytometric analysis of HepG2 cells using IL17RB antibody (bottom histogram) compared to a negative control (top histogram). FITC-conjugated goat-anti-rabbit secondary Ab was used for the analysis.



IHC analysis of FFPE human colon carcinoma stained with IL17RB antibody



Western blot analysis of IL17RB antibody and MDA-MB468 lysate.

## Handling & Storage

Handling	Aliquot the IL17RB antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Interleukin-17 receptor B
Application	ELISA, Flow Cytometry, IHC, WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human IL-17RB (aa 207-234).
Isotype	Ig
Purity Detail	Purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin sections) (1:10-1:50)Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q9NRM6
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 949 949 1129  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto: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)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)