

# Bcl-rambo (human) (CT) monoclonal antibody (Rocky-2)

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

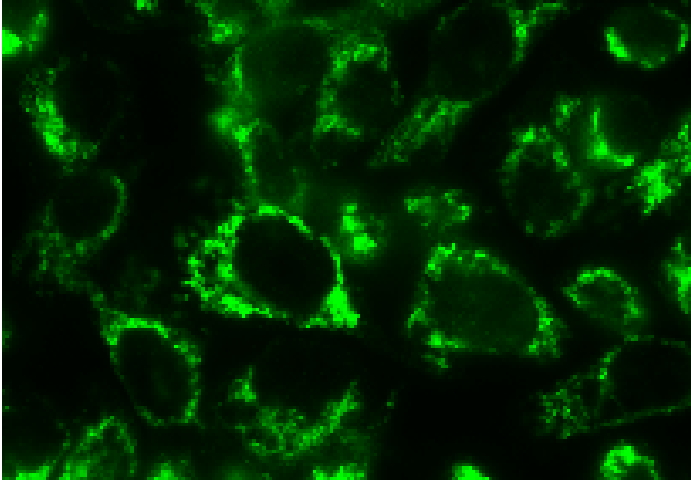
## Ordering Information

[Order Online »](#)

ALX-804-821-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Immunofluorescence microscopy showing the detection of endogenous Bcl-rambo in HELA cells with the MAb to Bcl-rambo (human) (CT) (Rocky-2) (Prod. No. ALX-804-821). **Method:** Following fixation with 2% paraformaldehyde/PBS, HELA cells were incubated with MAb to Bcl-rambo (human) (CT) (Rocky-2) (Prod. No. ALX-804-821) at a 1:300 dilution. An anti-mouse IgG antibody conjugated with Alexa Fluor 488 was used for detection.

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Bcl-2-like protein 13, Mil1 protein
Application	ICC, IP, WB
Application Notes	Detects endogenous protein by ICC, IP and WB.
Clone	Rocky-2
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human Bcl-rambo.
Isotype	IgG1
Recommendation Dilutions/Conditions	Immunocytochemistry (1:300)Immunoprecipitation (1:100)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human
Specificity	Recognizes the C-terminus of human Bcl-rambo (aa 224-459).
UniProt ID	Q9BXK5

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)