

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

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

Ordering Information

[Order Online »](#)

ALX-804-8210

100µg

Manuals, SDS & CofA

[View Online »](#)

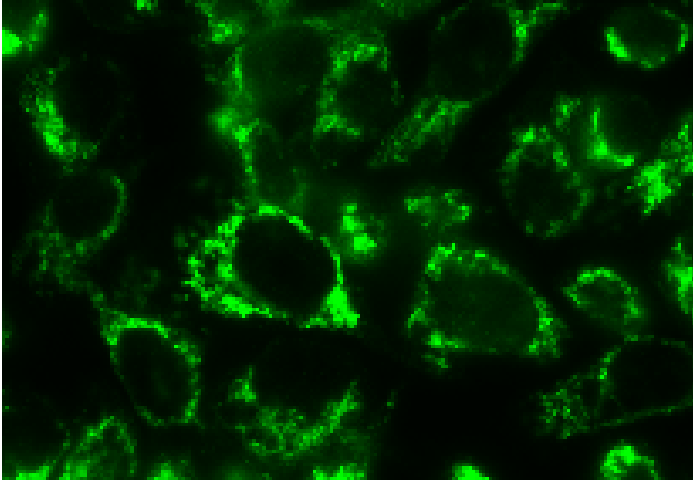


Figure: Immunofluorescence microscopy showing the detection of endogenous Bcl-rambo in HELA cells with the MAb to Bcl-rambo (human) (CT) (Rocky-2) (Prod. No. ALX-804-821). **Method:** Following fixation with 2% paraformaldehyde/PBS, HELA cells were incubated with MAb to Bcl-rambo (human) (CT) (Rocky-2) (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-2 (aa 224-459).

UniProt ID

Q9BXK5

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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