## 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 »** 

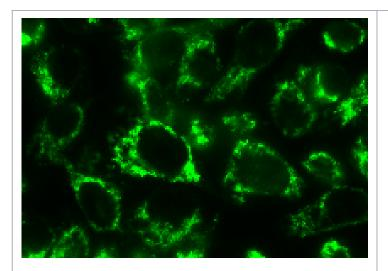


Figure: Immunofluorescence microscopy showing the detection of endogenous Bcl-rambo in HELA cells with the MAb to Bcl-rambo (human) (CT) (Rocky-2) (Prod. No. ALX-804-821). Method: Following fixation with 2% paraformal dehyde/PBS, HELA cells were incubated with MAb to Bcl-rambo (human) (CT) (Rocky-2)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

lgG1

**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

Last modified: May 29, 2024



info-