

# c-IAP2 monoclonal antibody (human) (16E-6-3)

The inhibitor of apoptosis protein (IAP) family consists of an evolutionarily conserved group of apoptosis inhibitors containing a conserved 70 aa BIR (baculovirus inhibitor repeat) domain. Human members of the family include c-IAP1, c-IAP2, XIAP, survivin, livin and NAIP. Overexpression of IAP family members, particularly survivin and livin, in cancer cell lines and primary tumors suggests an important role for these proteins in cancer progression. In general, the IAP proteins function through direct interactions to inhibit the activity of several caspases, including caspase-3, caspase-7 and caspase-9. In addition, binding of IAP family members to the mitochondrial protein Smac blocks its interaction with caspase-9, thereby allowing the processing and activation of the caspase.

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

Citations: 18

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-803-341-C100

100µg

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	HIAP-1, BIRC3, Baculoviral IAP repeat-containing protein-3, Cellular inhibitor of apoptosis-2, Human inhibitor of apoptosis protein-1, RNF49, API2
Application	WB
Clone	16E-6-3
Crossreactivity	Does not cross-react with mouse c-IAP2. Does not cross-react with human or mouse c-IAP1.
Formulation	Liquid. 0.2µm-filtered solution in PBS.
Host	Rat
Immunogen	GST-fusion product of the CARD domain of human c-IAP2.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q13489

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](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)