

# XIAP monoclonal antibody (2F1)

XIAP (X-linked inhibitor of apoptosis) belongs to the IAP (inhibitors of apoptosis) family, which is characterized by one or more BIR (baculovirus IAP repeat) domains. These proteins have multiple biological functions that include binding and inhibiting caspases, regulating cell cycle progression, and modulating receptor-mediated signal transduction. XIAP functions by direct binding and inhibition of several caspases, including caspase-3, caspase-7 and caspase-9.

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

Citations: 10 [View Online »](#)

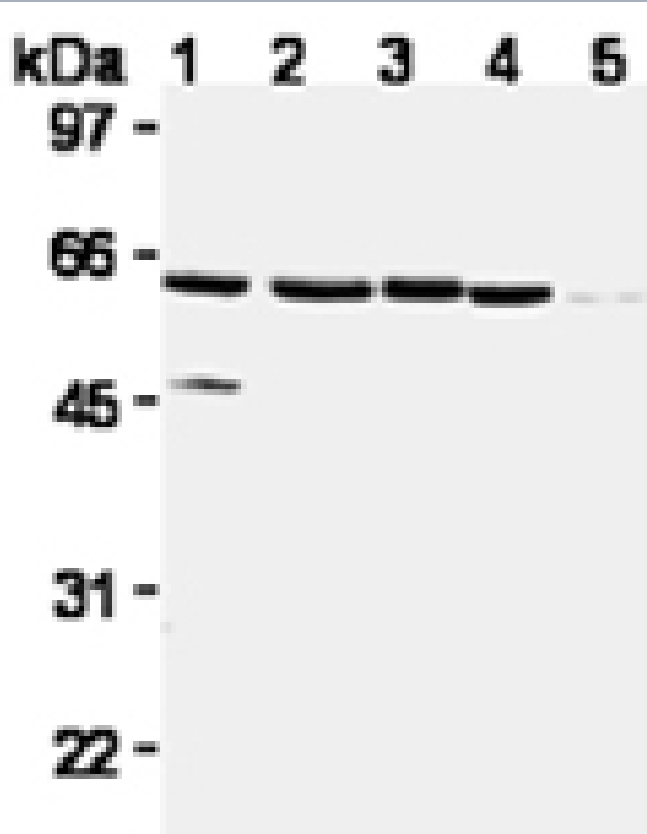
Ordering Information [Order Online »](#)

ADI-AAM-050-E	100µg
---------------	-------

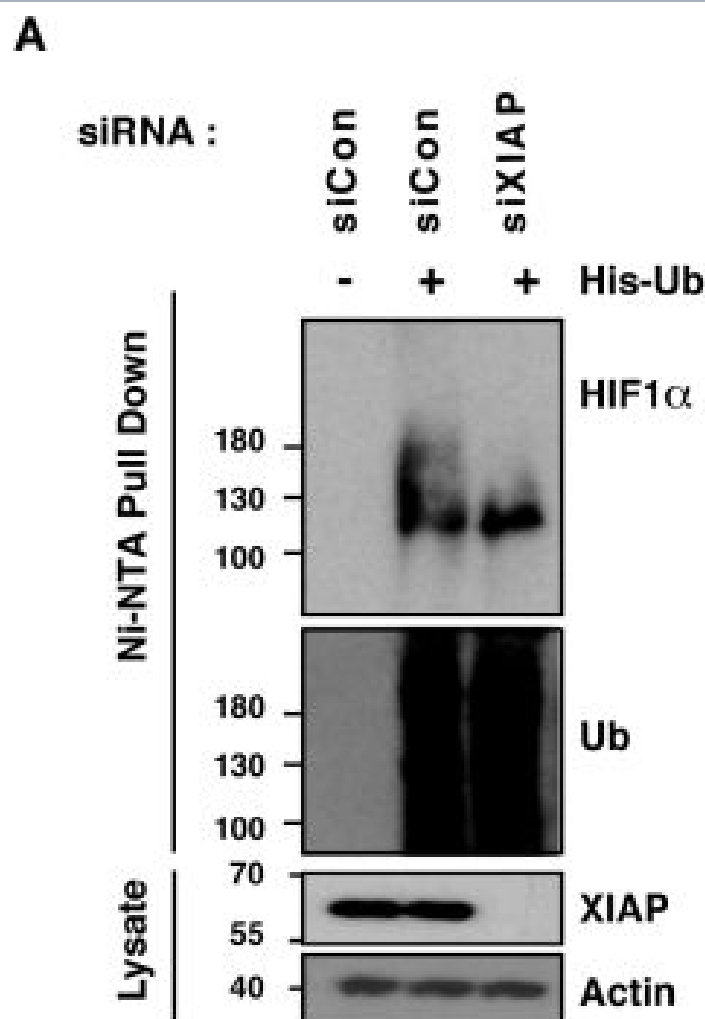
Manuals, SDS & CofA [View Online »](#)







Western blot analysis of Jurkat (1), Raji (2), HeLa (3), WR19L (4) and PC12 (5) cells, probed with XIAP mAb (2F1).



XIAP ubiquitinates HIF1α. (A) U2OS cells were transfected with siRNA oligonucleotides and His-tagged ubiquitin as indicated. Lysates were prepared and precipitated using nickel-conjugated beads followed by western blotting using the indicated antibodies. (B) HA-tagged plasmid encoding HIF1α was coexpressed in HEK293 cells with His-tagged ubiquitin (Ub), FLAG-XIAP (X) or FLAG-VHL (V) as indicated. Lysates were prepared 48 h post-transfection and ubiquitinated material was recovered by incubation with nickel-conjugated beads and analysed by western blotting. (C) Biotinylated-HIF1α was coexpressed in HEK293 cells with FLAG-tagged XIAP. HIF1α was recovered from lysates by incubation with streptavidin-coated beads and analysed using antibodies directed against specific ubiquitin chains. (D) U2OS cells were pre-treated with 2μM NSC697923 for 30 min before being exposed to 1% O<sub>2</sub> for 3 h. Cells were lysed in the presence of UBQ-TUBEs. Ubiquitin conjugates were purified using NiNTA agarose beads and analysed by immunoblotting using the indicated antibodies. (E) U2OS cells were pre-treated with 2 μM NSC697923 for 30 min before being exposed to 1% O<sub>2</sub> for 3 h. Cells were lysed in the presence of TAB2-TUBEs or TAB2-AA-TUBEs as indicated. Proteins conjugated to Lys63-linked ubiquitin chains were purified using Ni-NTA agarose beads and analysed by immunoblot.



## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      X-chromosome-linked inhibitor of apoptosis protein

Application      IHC, IP, WB

Application Notes      Detects a band of ~55kDa by Western blot.

Clone      2F1

Formulation      Liquid. In PBS, pH 7.2, containing 50% glycerol.

GenBank ID      AAC50373

Host      Mouse

Immunogen      Recombinant human XIAP.

Isotype      IgG1

Purity Detail      Protein A affinity purified.

**Recommendation**  
**Dilutions/Conditions**      Western Blot (1µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity      Human, Mouse

UniProt ID      P98170

**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)