

Livin/ML-IAP antibody (mAb)

Catalog No: 40958

RRID: AB_2793458

Clone: 88C570

Application(s): IHC, WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 31 kDa

Background: Livin (ML-IAP or BIRC7) protects against apoptosis induced by TNF or by chemical agents such as adriamycin, etoposide or staurosporine. Suppression of apoptosis is mediated by activation of MAPK8/JNK1, and possibly also MAPK9/JNK2. This activation depends on TAB1 and NR2C2/TAK1. *In vitro*, Livin (ML-IAP or BIRC7) inhibits Caspase-3 and proteolytic activation of pro-Caspase-9. Isoform 1 of Livin (ML-IAP or BIRC7) blocks staurosporine-induced apoptosis, while isoform 2 blocks etoposide-induced apoptosis.

Immunogen: This Livin/ML-IAP antibody was raised against a GST-fusion protein corresponding to human Livin/ML-IAP.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

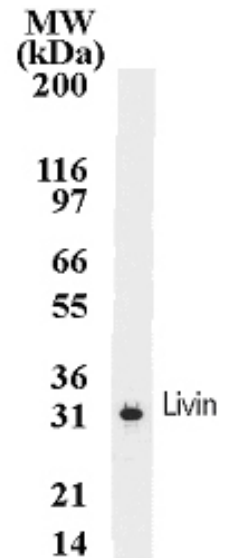
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Livin/ML-IAP mAb tested by Western blot. Livin detection by Western blot. The detection of Livin was performed using extract from a cell that had been transfected with human Livin cDNA. A protein of 31 kDa was detected using Livin/ML-IAP mAb at a 2 µg/ml dilution.