

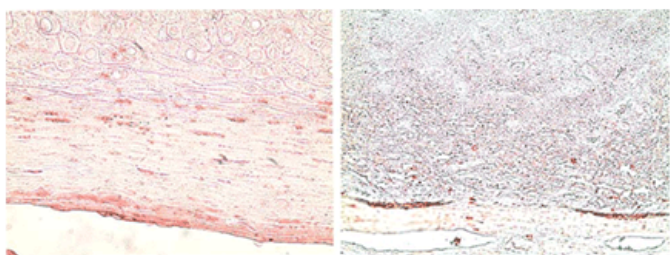
Survivin monoclonal antibody (60.11)

Survivin is a 16.5 kDa protein and the smallest inhibitor of apoptosis (IAP) so far identified. It is involved in the inhibition of apoptosis and cell division. Survivin expression has been reported at high levels in embryonic tissues, but at low or non-detectable levels in normal tissue. Survivin regulates the G2/M phase of the cell cycle by associating with the mitotic spindle microtubules and directly inhibits Caspase-3 and Caspase-7. Survivin is selectively expressed in the most common human cancers and is associated with clinical tumor progression. It has been proposed as a tumor marker for breast cancer, and Survivin expression has been correlated to clinical outcome in melanoma patients. Down-regulation or loss of Survivin is thought to inhibit the growth of tumor cells. Further, it has been indicated that Survivin epitopes may serve as important targets for anticancer immunotherapy approaches, and that Survivin is a rational target for apoptosis-based cancer therapy. It has also been proposed that Survivin may be used as a universal tumor antigen for immunotherapy.

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

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human tonsil tissue stained with Survivin monoclonal antibody (60.11), (Prod. No. ENZ-ABS715), at a 20 μ g/ml.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BIRC5, EPR1

Application IHC, WB

Clone 60.11

Formulation Liquid. In PBS containing 0.1% sodium azide.

Host Mouse

Immunogen Recombinant human Survivin protein (full length).

Isotype IgG2a

Purity Detail Protein G affinity purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rat

UniProt ID O15392

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)