

Smac/DIABLO

monoclonal antibody

(79-1-83)

Smac/DIABLO is one of the pro-apoptotic proteins normally found in the intermembrane space of the mitochondria. In response to apoptotic stimuli, Smac/DIABLO is released into the cytosol resulting in permeabilization of the mitochondrial membrane. Cytosolic Smac/DIABLO activates caspase-9 by binding IAP proteins. Changes in the expression level of Smac/DIABLO have been demonstrated in a number of human cancers including melanoma, ovary, and prostate.

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

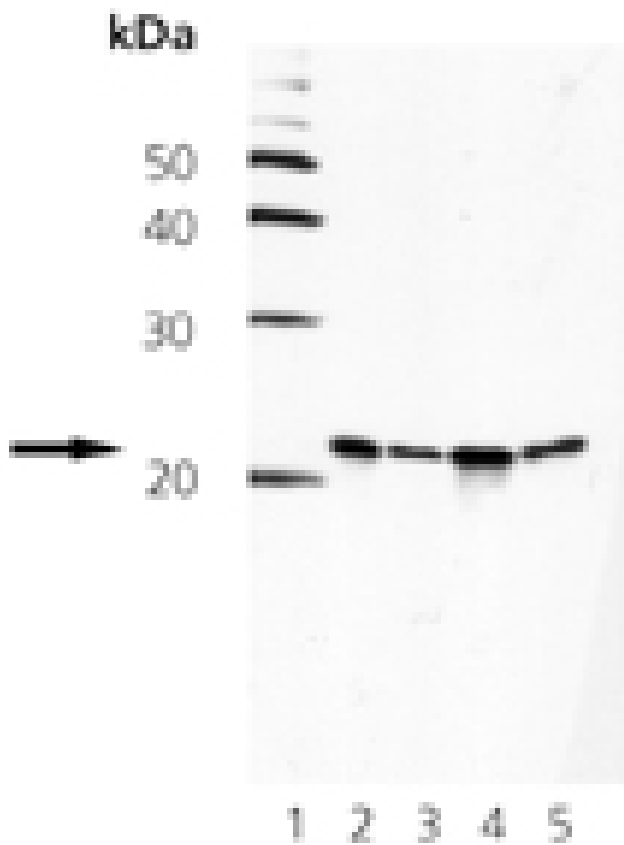
Ordering Information

[Order Online »](#)

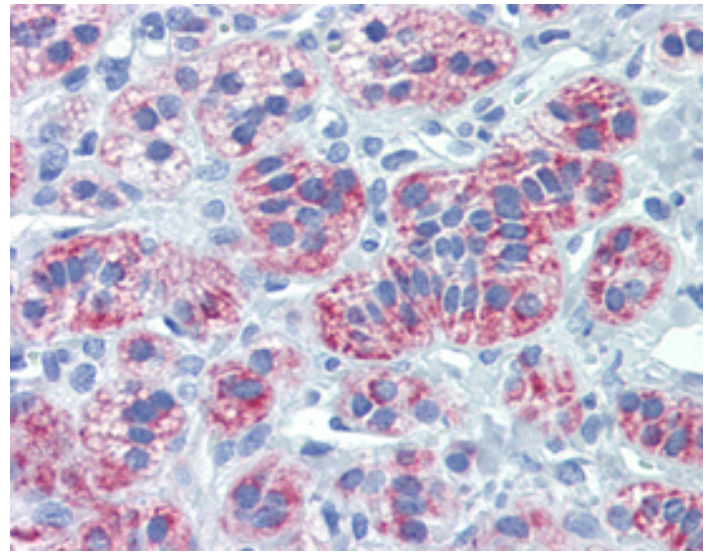
| | |
|-----------------|-------|
| ADI-905-684-100 | 100µg |
|-----------------|-------|

Manuals, SDS & CofA

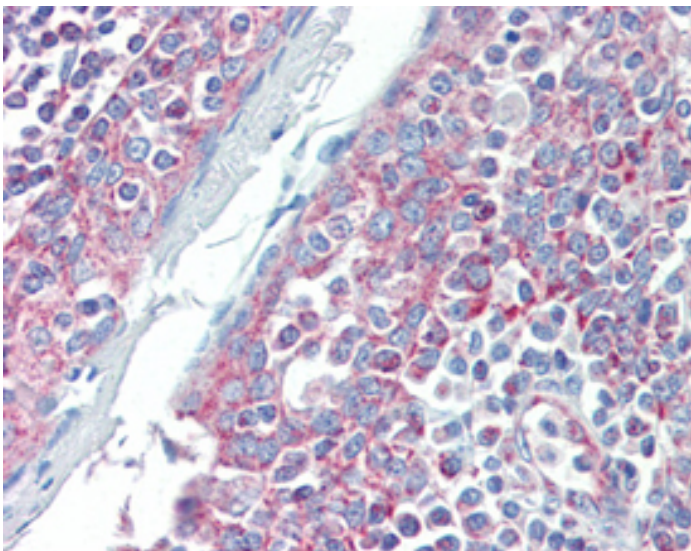
[View Online »](#)



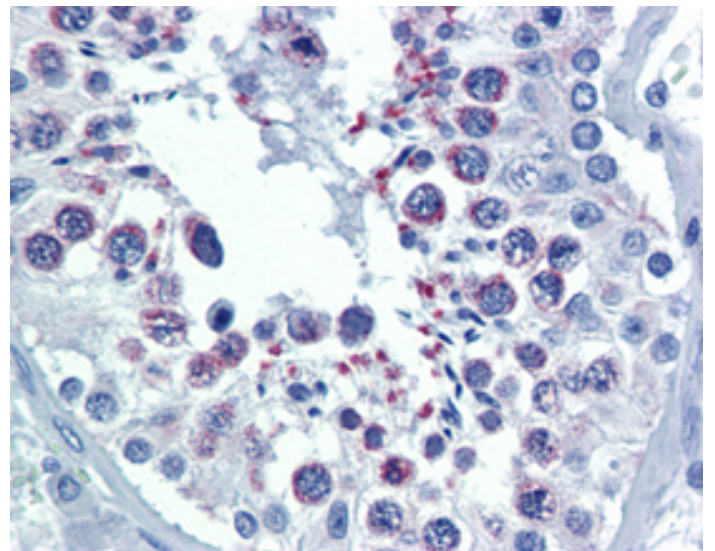
Western blot analysis of Smac/Diablo. Lane 1: MW marker, Lane 2: 293, Lane 3: HeLa, Lane 4: Jurkat, Lane 5: Vero.



IHC Analysis: Human adrenal tissue stained with Smac/DIABLO, mAb (79-1-83) at 10 µg/ml.



IHC Analysis: Human tonsil tissue stained with Smac/DIABLO, mAb (79-1-83) at 10 µg/ml.



IHC Analysis: Human testis tissue stained with Smac/DIABLO, mAb (79-1-83) at 10 µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Second mitochondria-derived activator of caspase

Application IHC (PS), WB

Clone 79-1-83

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.

Host Mouse

Immunogen Recombinant full length human Smac/DIABLO.

Isotype IgG2b

Purity Detail Protein G affinity purified.

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

Source Purified from mouse ascites.

Species Reactivity Human, Monkey

UniProt ID Q9NR28

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-
fr@enzolifesciences.com](mailto:info-fr@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)