

Caspase-8

monoclonal antibody

(5F7)

Caspase-8/FADD-like ICE (FLICE), is an ~55 kDa cytosolic protein belonging to the Fas/APO-1/CD95 death-inducing signaling complex. Although researchers have defined eight different isoforms of caspase-8 at the mRNA level (8/a-h), significant expression of only Caspase-8/a and /b occurs at the protein level in a variety of cell lines. Caspase-8 contains two death effector domains (DEDs) and exhibits sequence homology to both FADD, a signal transducer of Fas-induced apoptosis, and to the ICE/CED-3 family of cysteine proteases. Caspase-8 binds to FADD through its amino terminal DEDs and induces apoptosis when overexpressed. Activation of Caspase-8 involves a two-step proteolysis whereby the cleavage of Caspase-8 generates a 43 kDa fragment (p43) and a 12 kDa fragment further processed to 10 kDa, then the cleavage of receptor-associated p43 yields p26 and the release of the active site containing p18. A number of cytosolic factors inhibit the FADD-Caspase-8 containing signaling complex, including the ICE/CED-3 family inhibitor CrmA which inhibits Caspase-8 protease activity, and the short and long forms of FLIP that bind to the FADD-Caspase-8 complex.

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

Citations: 32

[View Online »](#)

Ordering Information

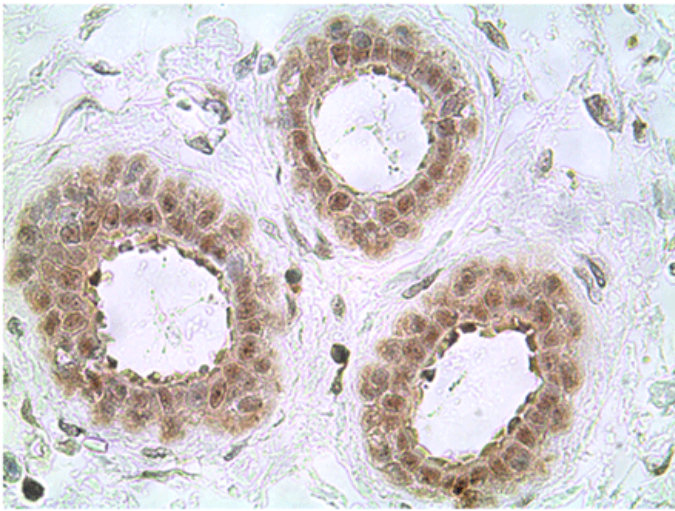
[Order Online »](#)

ADI-AAM-118-E	100µg
---------------	-------

Manuals, SDS & CofA

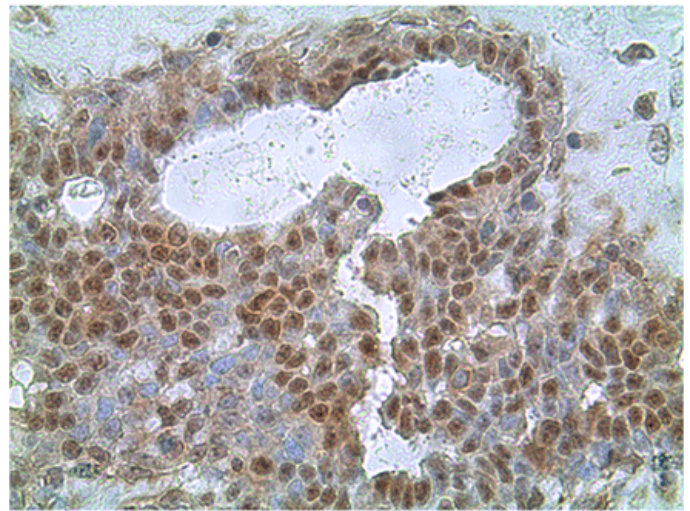
[View Online »](#)

63 x

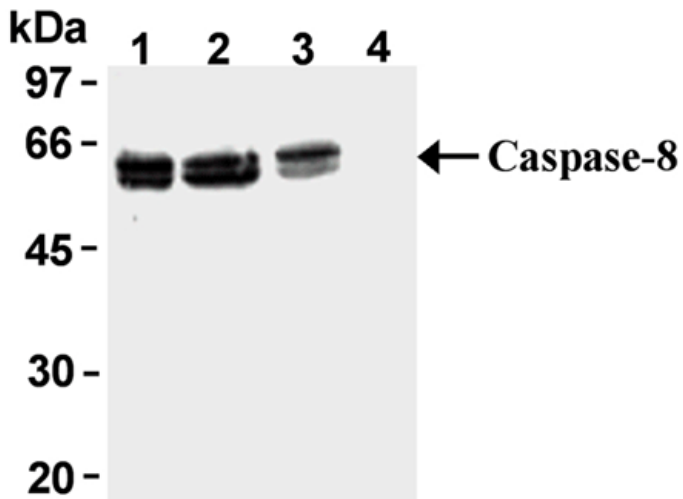


Immunohistochemistry staining of human breast tissue using anti-Caspase-8 antibody (prod.no. ADI-AAM-118 clone 5F7, 1:50) showed positive staining in both cytoplasm and nucleus of breast epithelial and myoepithelial cells. Image taken at magnification of 63x.

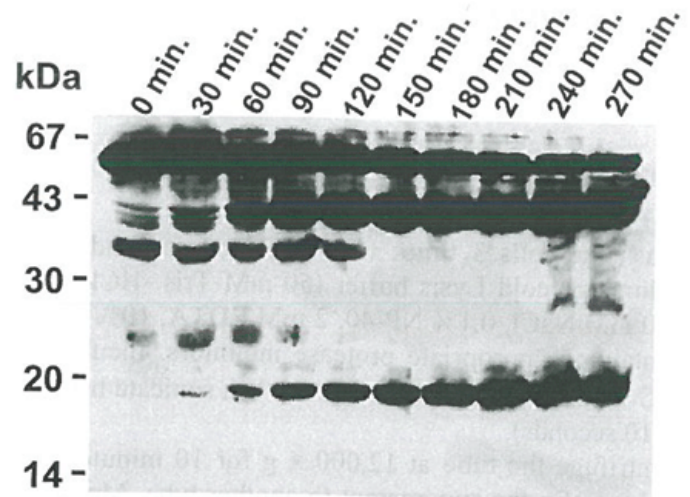
63 x



Immunohistochemistry staining of human breast tissue using anti-Caspase-8 antibody (prod.no. ADI-AAM-118 clone 5F7, 1:50) showed positive staining in both cytoplasm and nucleus of breast epithelial and myoepithelial cells. Image taken at magnification of 63x.



Western blot analysis of Jurkat (1), Raji (2), HeLa (3) and WR19L (4) probed with Caspase-8, mAb (5F7).



Western Blot analysis of Caspase-8 fragments in apoptosis induced Jurkat cells using Caspase-8, mAb (5F7) (Prod. No. ADI-AAM-118). The Jurkat cells were treated with a FAS monoclonal antibody for the indicated time

Long Term Storage	-20°C
Shipping	Blue Ice

Alternative Name	FLICE
Application	IHC (PS), IP, WB
Application Notes	Detects a band of ~55-57kDa by Western blot. Also detects the cleaved intermediate form (p41/43) and the activated large caspase-8 subunit p18 after apoptosis induction.
Clone	5F7
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Mouse
Immunogen	C-terminal fragment of recombinant human caspase-8 (amino acids 176-460).
Isotype	IgG2b
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:50)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q14790
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)