

c-FLIP polyclonal antibody

The cellular FLICE-like inhibitory protein, cFLIP, is a cellular homologue of the viral protein, vFLIP. cFLIP contains two death effector domains (DEDs) and is expressed as either a short or long splice variant called cFLIPs and cFLIPL, respectively. cFLIP acts as a caspase-8 dominant negative regulator, blocking apoptosis induced by the oligimerization of the adapter protein, FADD. cFLIP interacts with FADD and caspase-8 through their respective DEDs.

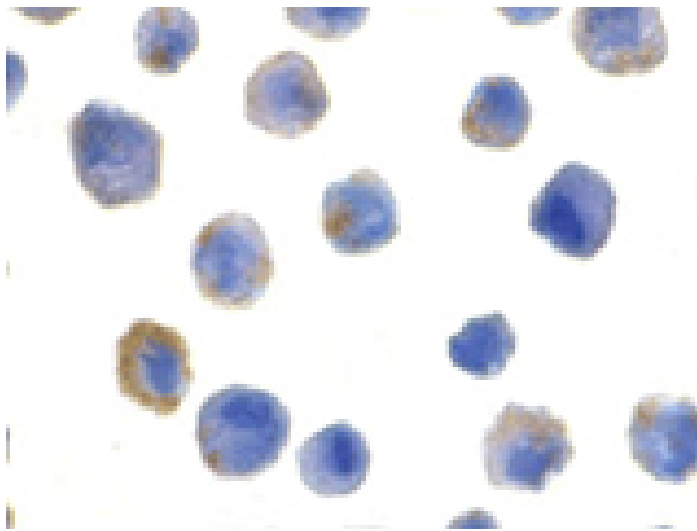
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 13 [View Online »](#)

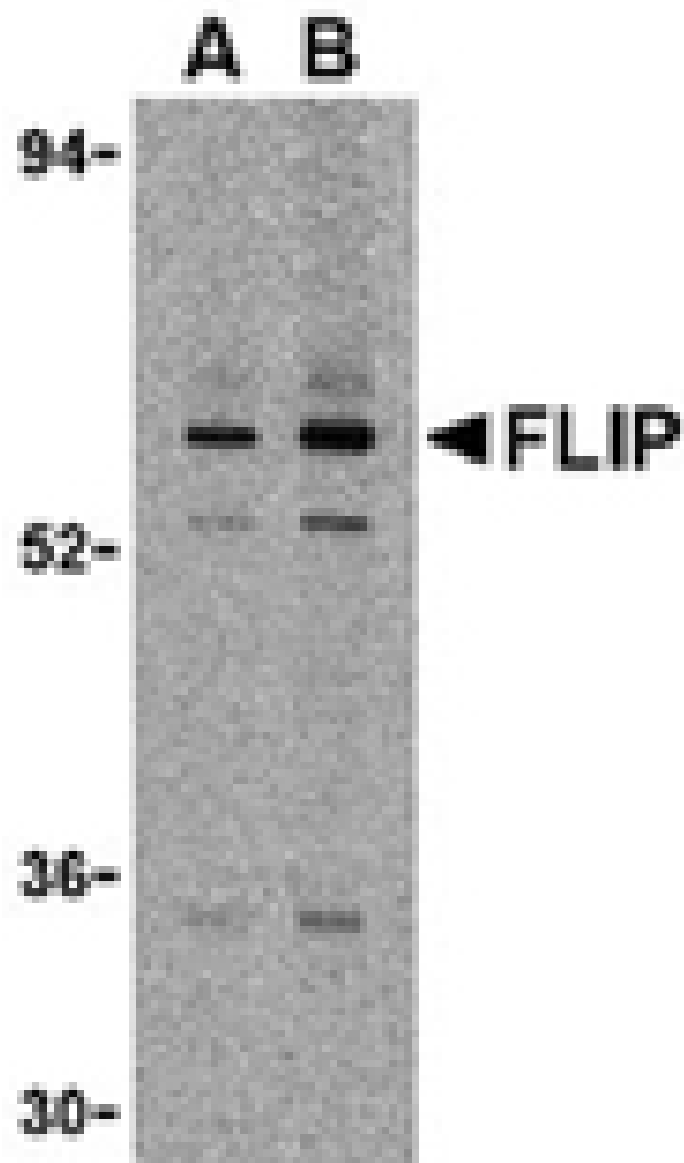
Ordering Information [Order Online »](#)

ADI-AAP-440-E	100µg
---------------	-------

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



Immunocytochemistry analysis of FLIP in HeLa cells with FLIP antibody at 5µg/ml.



Western blot analysis of FLIP in K562 cell lysate with FLIP antibody at (A) 1 and (B) 2µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CASPER, CLARP, I-FLICE, Usurpin

Application ICC, WB

Application Notes Detects a band of ~28-58kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

GenBank ID U97074, U97075

Host Rabbit

Immunogen FLIP antibody was raised against a peptide corresponding to 17 amino acids near the amino terminus of human FLIP. The sequence is identical in all FLIP splice variants. The immunogen is located within the first 50 amino acids of FLIP.

Purity Detail Peptide affinity purified.

Recommendation Immunocytochemistry (5µg/ml)Western Blot (0.5-1.0µg/ml)
Dilutions/Conditions Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID O15519

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)