

FLIP (human) monoclonal antibody (7F10)

FLIP contains two death effector domains (DEDs) and a caspase-like domain. FLIP interacts with the adapter protein, FADD, and caspase-8 and -10, and potentially inhibits apoptosis induced by death receptors CD95, DR3, TRAIL-R and TNFR1.

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

Citations: 57

[View Online »](#)

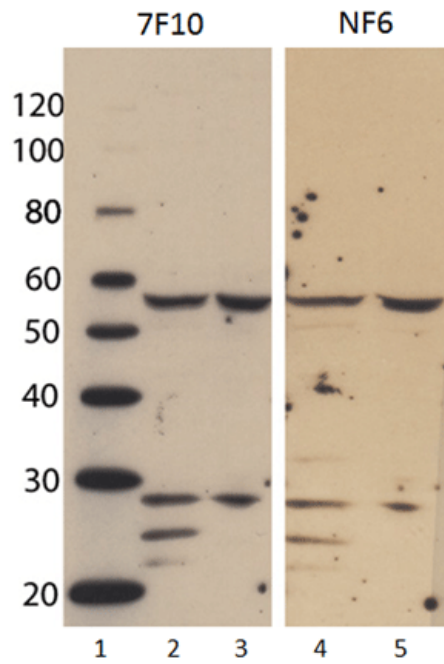
Ordering Information

[Order Online »](#)

ALX-804-961-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of FLIP (human), mAb (7F10) (Prod. No. ALX-804-961): Lane 1: MW marker, Lane 2&4: HeLa cell lysate, Lane 3&5: Jurkat cell lysate. The two membranes were blotted with 7F10 or NF6 clones at 2 μ g/ml separately. The result showed they detected very similar pattern in these cell lysates.

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	CASPER, CLARP, I-FLICE, Usurpin, CFLAR
Application	ELISA, Flow Cytometry, IHC, WB
Clone	7F10
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	Recombinant human FLIP.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Human
Technical Info / Product Notes	Cited samples: For an overview on cited samples please click here.
UniProt ID	O15519

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)