

FADD monoclonal antibody (1F7)

Fas-associating death domain containing protein (FADD)/MORT1 serves as a signal transducer of Fas-induced apoptosis. Through its amino-terminal death effector domain (DED), FADD binds to cytosolic proteins such as TRADD, a cytosolic death domaincontaining protein that associates with the cytosolic portion of TNF receptor, or MACH/FLICE/caspase-8, a protease with sequence homology to ICE-CED-3 protease family, in order to transduce the death signal.

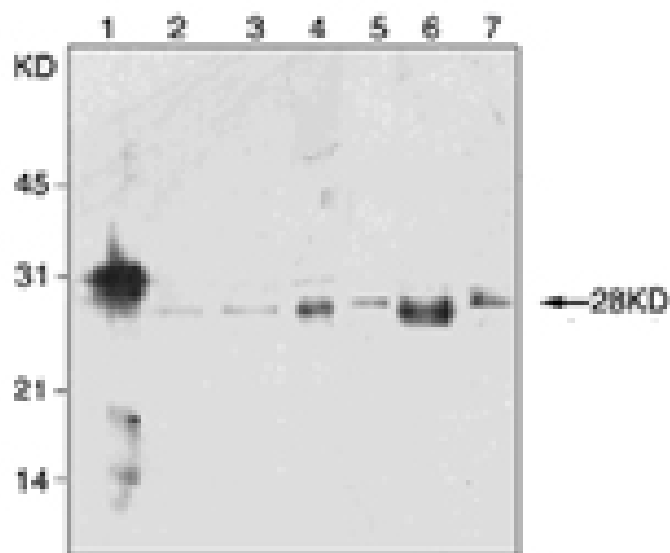
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 26 [View Online »](#)

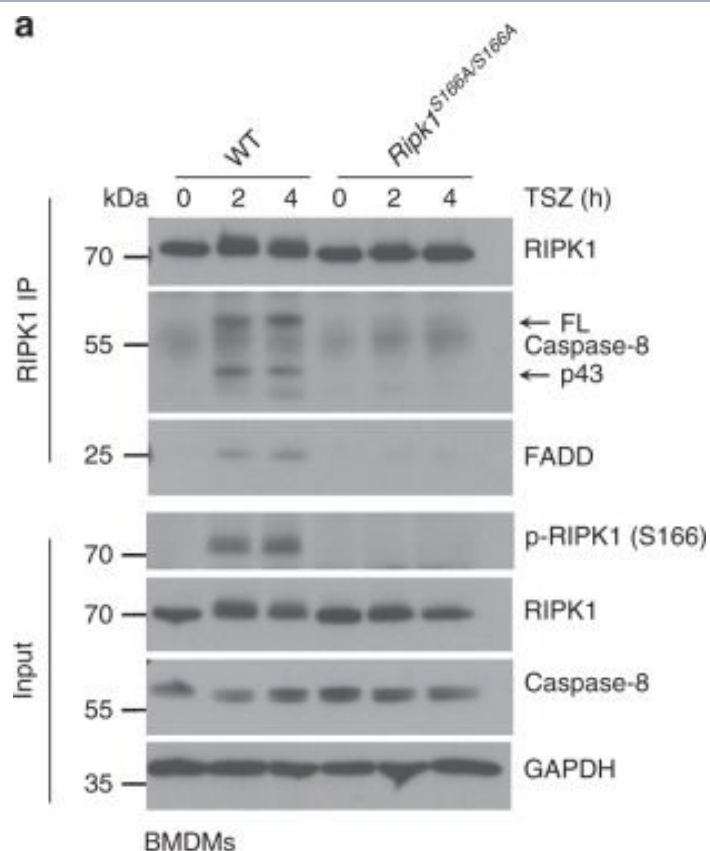
Ordering Information [Order Online »](#)

ADI-AAM-212-E	100µg
---------------	-------

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



Western blot probed with FADD mAb (1F7). Lane 1 FADD (recombinant) and cell lysates Jurkat (lane 2), Raji (lane 3), HeLa (lane 4), ZR75-1 (lane 5), A431 (lane 6), and HUC-Fm (lane 7).



S166 phosphorylation facilitates necrosome formation. a BMDMs from mice of the indicated genotypes were treated with TNF (T; 10 ng ml⁻¹), SMAC mimetic (S (Birinapant); 1 μM) and Z-VAD-FMK (Z) for 0, 2 and 4 h. RIPK1 immunoprecipitates and total cell lysates (input) were analyzed by immunoblotting with the indicated antibodies. b BMDMs from mice of the indicated genotypes were treated with TNF (T; 10 ng ml⁻¹), SMAC mimetic (S (Birinapant); 1 μM) and Z-VAD-FMK (Z) for 2 and 4 h. Lysates were processed under non-reducing conditions and analyzed for presence of MLKL by immunoblot. c BMDMs from mice of the indicated genotypes were treated with LPS (L; 100 ng ml⁻¹) and Z- VAD-FMK (Z) for 13 and 15 h. Lysates were processed under non-reducing or reducing conditions and analyzed by immunoblot using the indicated antibodies. d BMDMs from mice of the indicated genotypes were treated with Poly(I:C) (P; 0.5 μg ml⁻¹) and Z-VAD-FMK (Z) for 13 and 15 h. Lysates were processed under non-reducing or reducing conditions and analyzed by immunoblot using the indicated antibodies. a, c Representative data of three independent experiments is shown. b, d Representative data of two independent experiments is shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication:
Autophosphorylation at serine 166 regulates RIP kinase 1-mediated cell death and inflammation. *Nat Commun* (2020)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name MORT1, Fas (TNFRSF6)-associated via death domain

Application WB

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

Clone 1F7

Formulation Liquid. In PBS, pH 7.2, with 50% glycerol.

Host Mouse

Immunogen Recombinant human FADD.

Isotype IgG1

Purity Detail Protein A 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.

Species Reactivity Human, Mouse

UniProt ID Q13158

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



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com