

RAIDD (human) polyclonal antibody (AT100)

RAIDD (RIP-associated ICH-1 homologous protein with a death domain) is an adaptor molecule that mediates the action of cysteine proteases involved in apoptosis. In human cell lines, the caspase-2 adaptor RAIDD interacts selectively with caspase-2 through its caspase recruitment domain (CARD) and leads to caspase-2-dependent death. RAIDD overexpression induced caspase-2, CARD- and caspase-9-dependent apoptosis of PC12 cells and sympathetic neurons. Apoptosis correlated with the formation of discrete perinuclear aggregates. Both death and aggregates required the expression of full-length RAIDD. It has been reported that activation of caspase-2 occurs in a complex that contains the death domain-containing protein PIDD, whose expression is induced by p53, and the adaptor protein RAIDD.

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

Ordering Information

[Order Online »](#)

ALX-210-915-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

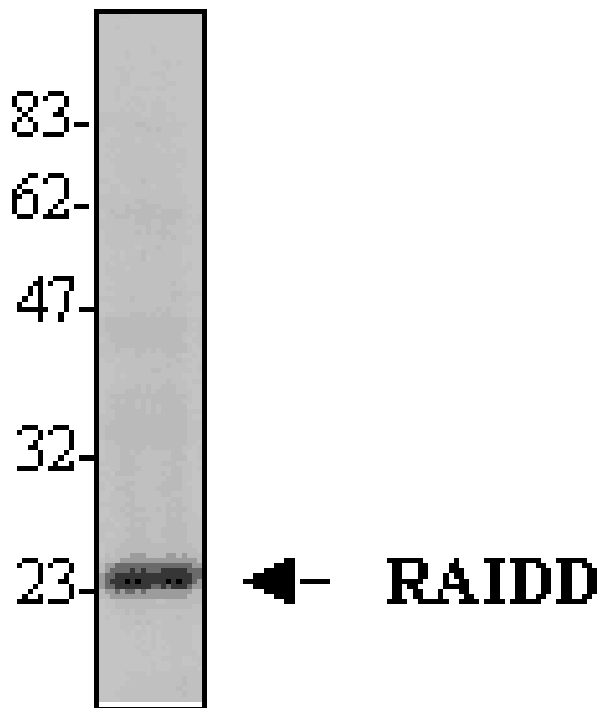


Figure: Detection of endogenous RAIDD (human) in HEK293T cells by the PAb to RAIDD (AT100) (Prod. No. ALX-210-915). **Method:** Total cell extracts (50µg) were run on a SDS-gel, and revealed with anti-RAIDD antibody (AT100, 1:1000, affinity purified antibody, in 5% milk powder) for 1 hour at RT followed by incubation with HRP-coupled anti-rabbit secondary antibody (1:4'000). The arrow indicates endogenous human RAIDD.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CRADD, Caspase and RIP adaptor with death domain, RIP-associated protein with a death domain
Application	IP, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human RAIDD (aa 1-199).
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	P78560

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
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-de@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