

# Viperin polyclonal antibody (AT131)

Viperin has been identified as a human cytomegalovirus and IFN- $\gamma$  inducible protein. Its expression is activated by both TLR3-dependent and RIG-I-dependent pathways. Viperin inhibits influenza A virus release of infected cells by altering plasma membrane fluidity and the formation of lipid rafts. It acts on the lipid rafts by decreasing the activity of the enzyme farnesyl diphosphate synthase (FPPS) which is essential for isoprenoid biosynthesis.

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

Citations: 4

[View Online »](#)

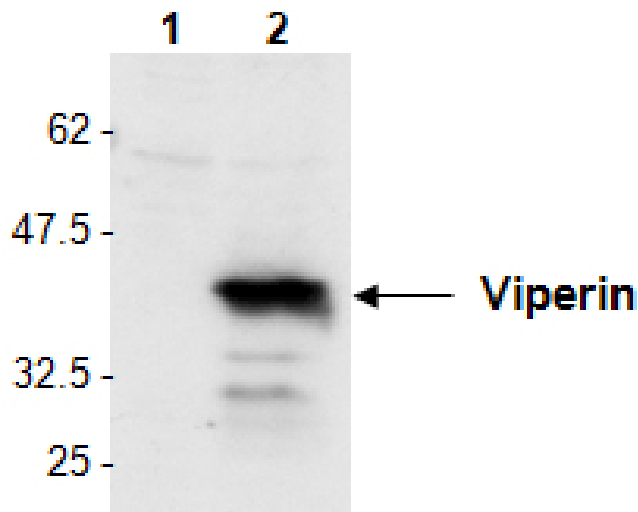
## Ordering Information

[Order Online »](#)

ALX-210-956-C100	100 $\mu$ g
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Detection of mouse Viperin in L929 cells by Western Blot using PAb to Viperin (AT131) (Prod. No.

[ALX-210-956](#)) **Method:** Cell extracts from L929 cell either unstimulated (lane 1) or stimulated for 24h with poly(dA.dT) poly(dT.dA) at 3µg/ml (lane 2) were resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and incubated with the PAb to Viperin (AT131) at a 1:1000 dilution for 1 hour. Proteins were visualized using a peroxidase-conjugated polyclonal antibody to rabbit IgG and a chemiluminescence detection system. """, Viperin polyclonal antibody (AT131) Western blot, Viperin polyclonal antibody (AT131) Western blot, yes ALX-215-001, 215-

001.gif, <http://static.enzolifesciences.com/fileadmin/files/image/215-001.gif>,

**Figure:** Peptide competition. Only the phosphopeptide corresponding to tau [pS<sup>262</sup>] blocks the antibody signal

# 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	Cig5, RSAD2, Cytomegalovirus-induced gene 5 protein, Radical S-adenosyl methionine domain-containing protein 2, Virus inhibitory protein, endoplasmic reticulum-associated, interferon-inducible
Application	IP, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa near the C-terminus of human viperin.
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:200)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, Mouse
UniProt ID	Q8WXG1
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)