

# Caveolin-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-241-C100	100µg
------------------	-------

Manuals, SDS & CofA

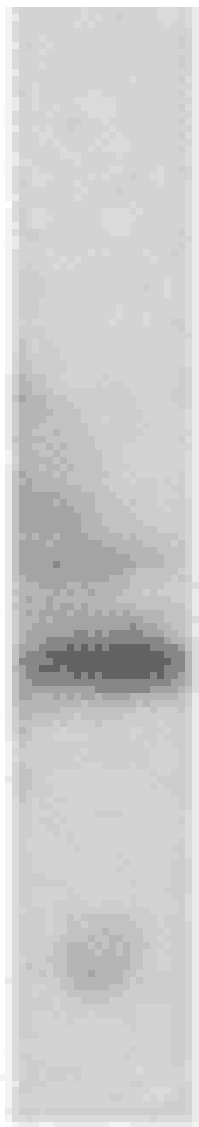
[View Online »](#)

38 kDa-

25 kDa-

20 kDa-

14 kDa-



**Figure:** Western blot of caveolin-3 on rat cardiac muscle protein extract using PAb to Caveolin-3 (Prod. No. ALX-210-241).

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** ICC, IF, IHC, IP, WB

**Application Notes** Detects a band of ~21kDa by Western blot.

**Crossreactivity** Does not cross-react with caveolin-1 or -2.

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 1-18 (M1MTEEHTDLEARIIKDIH18C) of rat caveolin-3.

**Purity Detail** Epitope affinity-purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:100-200)Immunofluorescence (1:20)Western Blot (1µg/ml)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** P51638

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



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 940 948 9120  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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