

Caveolin-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-2410	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

38 kDa-

25 kDa-

20 kDa-

14 kDa-

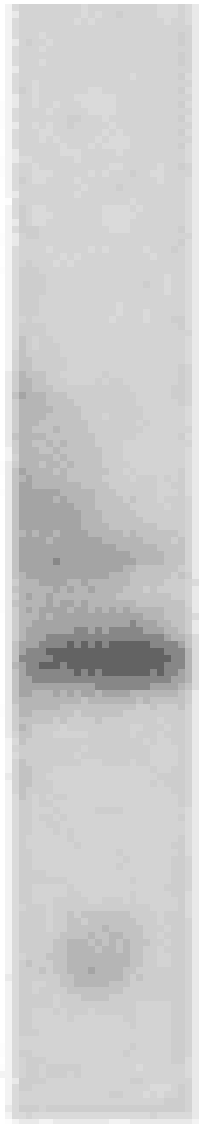


Figure: Western blot of caveolin-3 on rat cardiac muscle protein extract using PAb to Caveolin-3 (Prod. No. ALX-210-241).

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