

Actin polyclonal antibody

Actin is one of the major proteins (~42 kDa) of the cytoskeleton. It regulates contractile potential in muscle tissues and helps to control the shape as well as the motility of non-muscle cells. Actin is expressed in all eukaryotic cells which makes it a useful loading control in Western blotting experiments. In higher eukaryotes, there are six isoforms of actin: Three alpha-actins (α -skeletal, α -cardiac, α -smooth muscle), one β -actin (β -non-muscle) and two γ -actins (γ -smooth muscle and γ -non-muscle). Actin isoforms show >90% overall sequence homology.

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

Citations: 4

[View Online »](#)

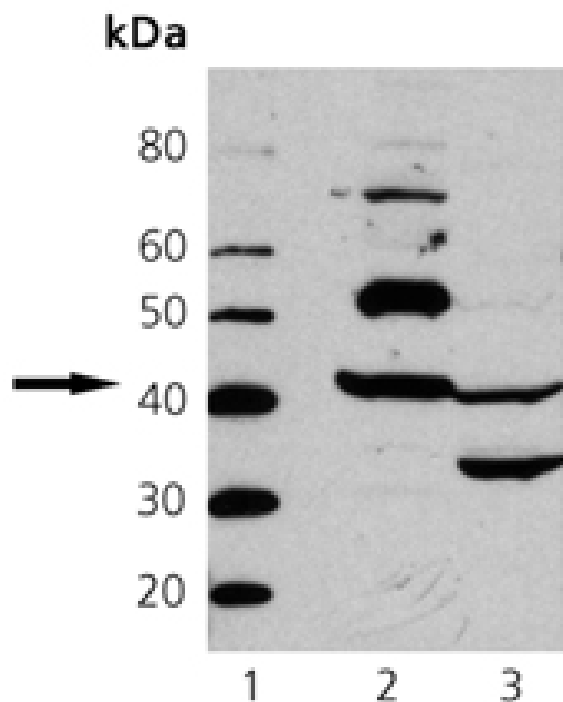
Ordering Information

[Order Online »](#)

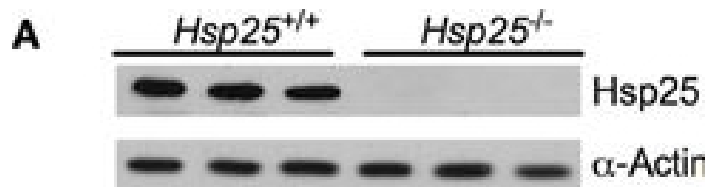
ADI-CSA-400-D	50 μ g
ADI-CSA-400-F	200 μ g

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Actin pAb: Lane 1: MWM, Lane 2: Mouse Heart Tissue Extract, Lane 3: Chicken Heart Tissue Extract.



(A) Gastrocnemius muscle HSP25 and actin protein levels in Hsp25^{+/+} and Hsp25^{-/-} mice verifying the absence of HSP25. (B) HSP25 protein levels in young and old Hsp25^{+/+} mice under sedentary conditions or following 14 days of running wheel activity.

#Significantly different from sedentary ($P < 0.0001$),
†significant main effect of age ($P < 0.0001$).

Representative western blot showing HSP25 in sedentary and running wheel young and old Hsp25^{+/+} mice. Data are presented as means \pm SE, $n = 8$ for all groups.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Effect of HSP25 loss on muscle contractile function and running wheel activity in young and old mice. *Front Physiol* (2014)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name ACTA, ACTSA

Application WB

Application Notes Detects a band of ~42kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.

Host Rabbit

Immunogen Bovine skeletal muscle Actin.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL) 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 Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID P68133

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

.