

p23 polyclonal antibody

p23 is a ubiquitous, highly conserved co-chaperone for Hsp90 that participates in the folding of a number of cell regulatory proteins including progesterone and glucocorticoid receptors, HSF1, and telomerase. p23 is thought to modulate Hsp90 activity in the last stages of the chaperoning pathway by binding and stabilizing Hsp90 in its ATP-bound state.

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

Citations: 7

[View Online »](#)

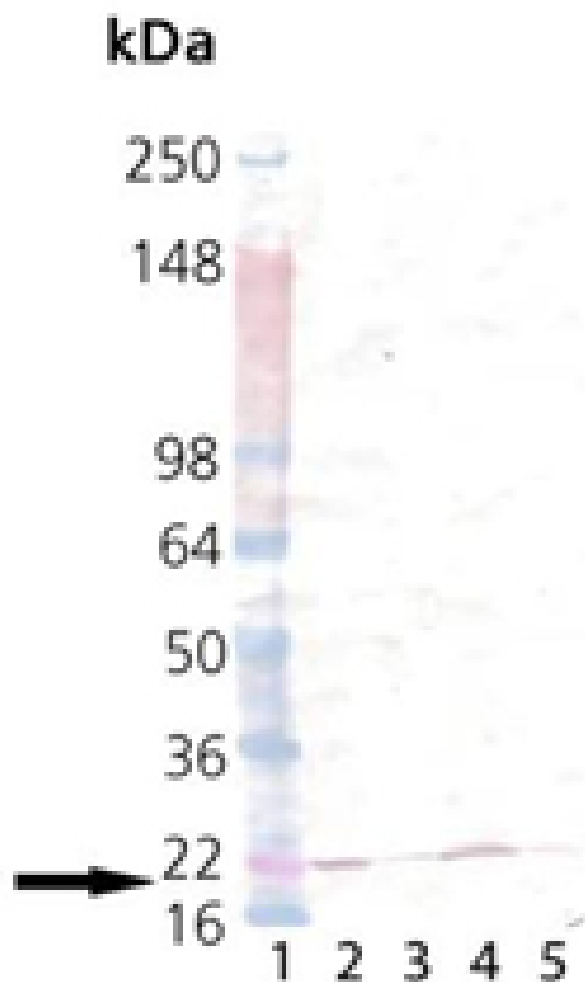
Ordering Information

[Order Online »](#)

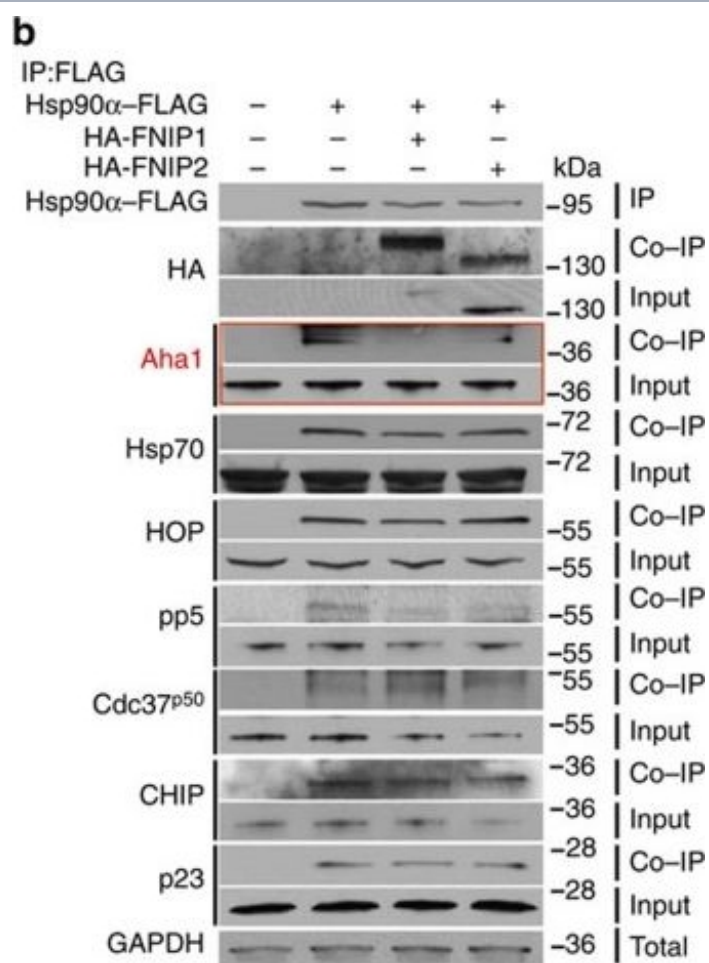
ADI-SPA-670-D	50µl
ADI-SPA-670-F	200µl

Manuals, SDS & CofA

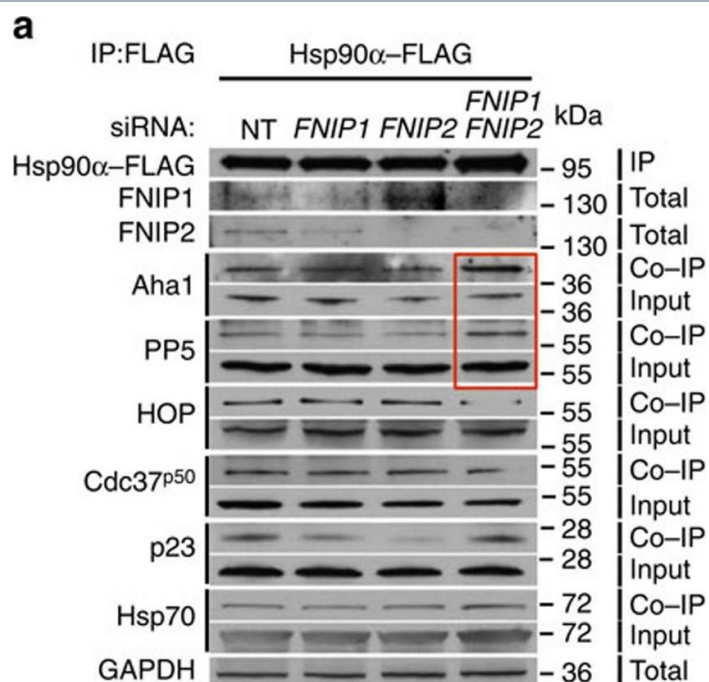
[View Online »](#)



Western blot analysis of p23: Lane 1: MW, Lane 2: p23 (Prod. No. ADI-SPP-670), Lane 3: HeLa (Heat Shocked), Lane 4: mouse testes, and Lane 5: rat testes probed with p23, pAb (Prod. No. ADI-SPA-670).



FNIPs compete with the Aha1 co-chaperone for binding to Hsp90. (a) Hsp90α-FLAG was transiently expressed in HEK293 cells for 24 h followed by siRNA knockdown of FNIP1 and/or FNIP2. Hsp90α-FLAG was immunoprecipitated (IP) and co-IP of the co-chaperones was assessed by immunoblotting. Densitometry of the western blotting for FNIPs is represented as mean±s.d. A Student's t-test was performed to assess statistical significance (*P<0.05, **P<0.005 and ***P<0.0005). (b) HEK293 cells were transiently co-transfected with Hsp90α-FLAG and HA-FNIP1 or HA-FNIP2. Hsp90α-FLAG was isolated and co-IP of co-chaperones examined by immunoblotting. (c) HA-FNIP1, HA-FNIP1-D and HA-FNIP2 inhibited Hsp90α-HA ATPase activity after 30 min. Addition of 1.3 μM Aha1-FLAG stimulated the ATPase activity. All the data represent mean±s.d. A Student's t-test was performed to assess statistical significance (**P<0.005 and ****P<0.0001). (d) FNIP1 and Aha1 compete for binding to Hsp90α. FNIP1-D-His6 was attached to Ni-NTA agarose and then incubated with Hsp90α. Ni-NTA agarose was then washed and incubated with the indicated amounts of Aha1-FLAG. (e) Aha1-FLAG attached to anti-FLAG M2 affinity gel was incubated with Hsp90α initially and then washed and incubated with indicated amounts of the FNIP1-D-His6.



FNIPs compete with the Aha1 co-chaperone for binding to Hsp90. (a) Hsp90α-FLAG was transiently expressed in HEK2993 cells for 24 h followed by siRNA knockdown of FNIP1 and/or FNIP2. Hsp90α-FLAG was immunoprecipitated (IP) and co-IP of the co-chaperones was assessed by immunoblotting. Densitometry of the western blotting for FNIPs is represented as mean±s.d. A Student's t-test was performed to assess statistical significance (*P<0.05, **P<0.005 and ***P<0.0005). (b) HEK2993 cells were transiently co-transfected with Hsp90α-FLAG and HA-FNIP1 or HA-FNIP2. Hsp90α-FLAG was isolated and co-IP of co-chaperones examined by immunoblotting. (c) HA-FNIP1, HA-FNIP1-D and HA-FNIP2 inhibited Hsp90α-HA ATPase activity after 30 min. Addition of 1.3 μM Aha1-FLAG stimulated the ATPase activity. All the data represent mean±s.d. A Student's t-test was performed to assess statistical significance (**P<0.005 and ****P<0.0001). (d) FNIP1 and Aha1 compete for binding to Hsp90α. FNIP1-D-His6 was attached to Ni-NTA agarose and then incubated with Hsp90α. Ni-NTA agarose was then washed and incubated with the indicated amounts of Aha1-FLAG. (e) Aha1-FLAG attached to anti-FLAG M2 affinity gel was incubated with Hsp90α initially and then washed and incubated with indicated amounts of the FNIP1-D-His6.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: The FNIP co-chaperones decelerate the Hsp90 chaperone cycle and enhance drug binding. *Nat Commun* (2016)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Prostaglandin E synthase 3, HSP90 Co-chaperone

Application WB

Application Notes Detects a band of ~23kDa by Western blot.

Formulation Serum.

Host Rabbit

Immunogen Recombinant human p23.

Purity Detail Antiserum.

Recommendation Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available
Dilutions/Conditions for all applications. Optimal conditions must be determined individually for each application.

Source Whole rabbit serum.

Species Reactivity Human, Mouse, Rat

UniProt ID Q15185

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



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)