

FKBP59 monoclonal antibody (KN382/EC1)

FKBP59, also known as Hsp56/p59 is an immunophilin that is a component of Hsp70/Hsp90/steroid receptor complexes. Expression of FKBP59 is increased in response to heat shock and ischemic stress, and is thought to play a role in the folding and cytoplasmic/nuclear distribution of steroid receptors.

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

Citations: 3

[View Online »](#)

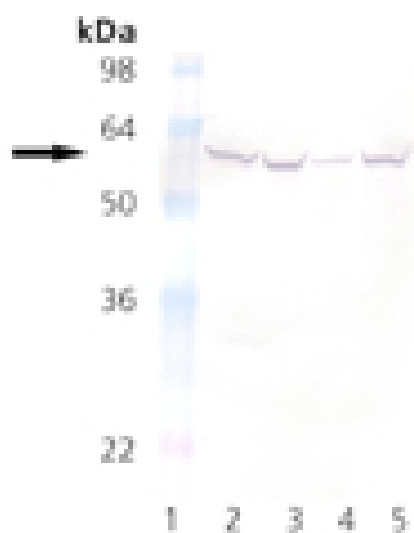
Ordering Information

[Order Online »](#)

ADI-SRA-1400-D	50µg
ADI-SRA-1400-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of FKBP59 mAb (KN382/EC1) (Prod. No. ADI-SRA-1400): Lane 1: MW marker, Lane 2: HeLa Cell Lysate(heat shocked), Lane 3: MDBK Cell Lysate (heat shocked), Lane 4: VERO Cell Lysate (heat shocked), Lane 5: RK-13 Cell Lysate (heat shocked).

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Product Details

Alternative Name	HSP56, Heat shock protein 56, FK506 binding protein 1A
Application	IHC, IP, WB
Application Notes	Detects a band of ~59kDa by Western blot.
Clone	KN382/EC1
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M84474
Host	Mouse
Immunogen	Partially purified 9S rabbit uterine progesterone receptor.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (10µg/ml)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from mouse ascites.
Species Reactivity	Bovine, Guinea pig, Hamster, Human, Mink, Monkey, Mouse, Rabbit

Technical Info / Product Notes

Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000004478).

UniProt ID

P27124

Worry-free Guarantee

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

Last modified: May 29, 2024



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)