

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 »](#)

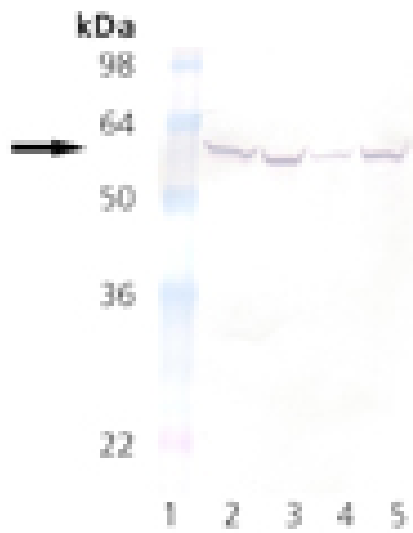
Ordering Information

[Order Online »](#)

ADI-SRA-1400-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 & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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