

FKBP12 polyclonal antibody

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-142-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

46 kDa

30 kDa

21 kDa

14 kDa



FKBP12

Figure: Western blot of purified FKBP12 after immunoprecipitation on a 15% gel using PAb to FKBP12 (Prod. No. [ALX-210-142](#)).^{""},FKBP12 polyclonal antibody Western blot,FKBP12 polyclonal antibody Western blot,yes
ALX-210-154,210-154.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-154.gif>,Figure 1 shows a Western blot of GRb on COS-74 whole cell extract using Prod. No. [ALX-210-154](#).^{""""""}

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 FK506 binding protein (12kDa)

Application IHC (FS), IHC (PS), IP, WB

Application Notes Detects a band of ~12kDa by Western blot.

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 1-13 (G1VQVETISPGDGR13) of human FKBP12 (FK506 binding protein 12kDa).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 2µg/ml)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 P62942

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
[usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
[fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
[fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
[de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

Last modified: May 29, 2024