

FKBP12 polyclonal antibody

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-1420	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](#)).¹,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-](http://static.enzolifesciences.com/fileadmin/files/image/210-154.gif)

154.gif,Figure 1 shows a Western blot of GRb on COS-74 whole cell extract using Prod. No. [ALX-210-154](#).²

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

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

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

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

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

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

Last modified: May 29, 2024