

PARP-2 (mouse) polyclonal antibody

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-8990	100µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for 1 week when stored at +4°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Poly(ADP-ribose) polymerase-2
Application	ICC, IHC, IP, WB
Crossreactivity	Weakly cross-reacts with human PARP-2.
Formulation	Liquid. Antiserum containing 0.1% sodium azide.
Host	Rabbit
Immunogen	Recombinant mouse PARP-2 (poly(ADP-ribose) polymerase-2) fragment (aa 1-207) produced in Sf21 cells.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:500-1:1,000)Immunohistochemistry (1:500-1:1,000)Immunoprecipitation (2µl per 100µg nuclear extract)Western Blot (1:2,000-1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. Immunohistochemistry (frozen sections, or PFA-fixed cryosections)
Species Reactivity	Mouse
UniProt ID	O88554
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)