

# 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

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Poly(ADP-ribose) polymerase-2
<b>Application</b>	ICC, IHC, IP, WB
<b>Crossreactivity</b>	Weakly cross-reacts with human PARP-2.
<b>Formulation</b>	Liquid. Antiserum containing 0.1% sodium azide.
<b>Host</b>	Rabbit
<b>Immunogen</b>	Recombinant mouse PARP-2 (poly(ADP-ribose) polymerase-2) fragment (aa 1-207) produced in Sf21 cells.
<b>Recommendation Dilutions/Conditions</b>	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)
<b>Species Reactivity</b>	Mouse
<b>UniProt ID</b>	O88554
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)