

PARP-2 monoclonal antibody (4G8)

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

Citations: 11

[View Online »](#)

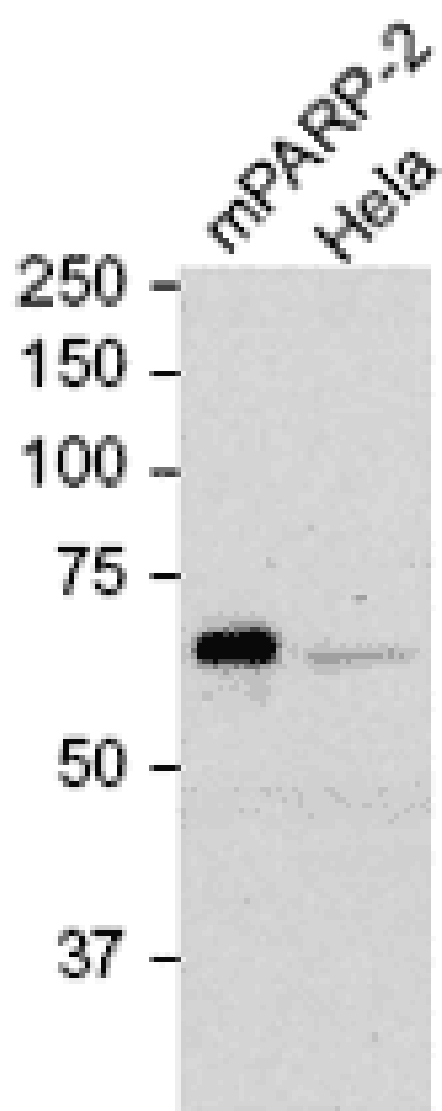
Ordering Information

[Order Online »](#)

ALX-804-639-L001	1ml
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Poly(ADP-ribose) polymerase-2
Application	ELISA, WB
Application Notes	Detects a band of ~62kDa by Western blot.
Clone	4G8
Formulation	Liquid. Tissue culture supernatant containing 0.1% sodium azide.
Host	Mouse
Immunogen	Recombinant mouse PARP-2 (poly(ADP-ribose) polymerase-2).
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (1:500)Western Blot (1:20-1:50)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
Specificity	Recognizes an epitope in domain E of PARP-2.
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

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