

PARP-10 (human) monoclonal antibody (5H11)

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

Citations: 1

[View Online »](#)

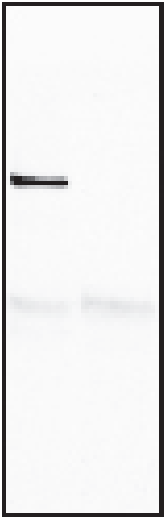
Ordering Information

[Order Online »](#)

ALX-804-626-L001	1ml
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

PARP10 →  170
130

pCMV - +
pCMV-PARP10 + -

Figure: Detection of PARP-10 using PARP-10 (human), mAb (5H11) (Prod. No. ALX-804-626) **Method:** HEK293 cells were transfected with a CMV-PARP10 or CMV vector (control). Frackelton extracts were used for Western blotting. The signal was detected using anti-rat-HRP and enhanced chemiluminescence.

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Poly(ADP-ribose) polymerase-10
Application	ICC, IP, WB
Clone	5H11
Formulation	Liquid. Tissue culture supernatant (TCS) containing 0.1% sodium azide.
Host	Rat
Immunogen	Recombinant human PARP-10 (poly(ADP-ribose) polymerase-10) (aa 1-907).
Isotype	IgG1
Recommendation Dilutions/Conditions	Immunocytochemistry (1:10)Immunoprecipitation (1:10)Western Blot (1:50)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q53GL7
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)