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

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

Ordering Information

ALX-804-626-L001

1ml

Manuals, SDS & CofA

View Online »

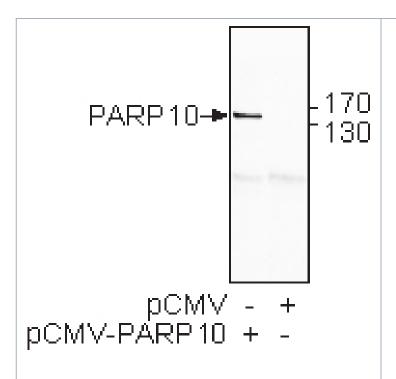


Figure: Detection of PARP-10 using PARP-10 (human), mAb (5H11) (Prod. No. ALX-804-626) **Method:** HEK293 cells were transfected with a CMV-PARP10 or CMV vector (control). Frackelton extracts were used for Western blotting. The signal was detected using anti-rat-HRP and enhanced chemiluminescence.

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).

lgG1

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com