

# Poly(ADP-ribose) monoclonal antibody

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

Citations: 23

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA216-0100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** PAR

**Application** ELISA, ICC, IP, WB

**Clone** 10HA

**Formulation** Liquid. In TBS.

**Host** Mouse

**Immunogen** Purified poly(ADP-ribose) polymer (chain length of 2-50 units).

**Isotype** IgG3

**Purity Detail** Purified from mouse ascites.

**Recommendation Dilutions/Conditions** Western Blot (1:1,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Species independent

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

Last modified: February 4, 2025



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)