

Poly(ADP-ribose), standard

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-202-043-C001	1µg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for one week when stored at +4°C. Stable for at least 2 years after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PAR
Application Notes	Immunoblotting standard for the detection of poly(ADP-ribose) (PAR) of 3-200 bases in size synthesized <i>in vitro</i> or <i>in vivo</i> .
Formulation	Liquid. In water.
MW	1'100 to 275'000 depending on the number of units of ADP-ribose.
Purity	≥95% (HPLC of ADP-ribose RNA and DNA measurement)
Quantity	1µg representing 1.8 nanomoles of 2-300 units of ADP-ribose.
Source	Released from automodified PARP-1 by base treatment and purified by boronate affinity.



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-
fr@enzolifesciences.com](mailto:info-fr@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)