

PARP-1 (human), (recombinant) (high purity)

Citations: 14

[View Online »](#)

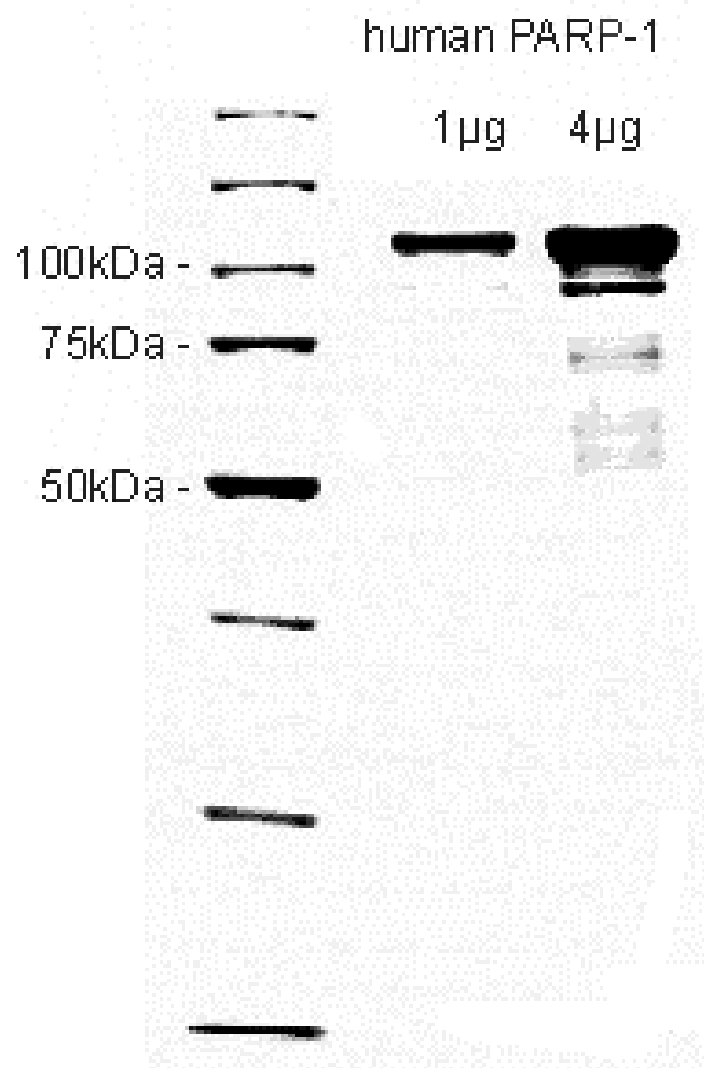
Ordering Information

[Order Online »](#)

ALX-201-063-C020	20µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Quality Control: 10% SDS-Polyacrylamide gel.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Poly(ADP-ribose) polymerase 1

Concentration 1mg/ml

Formulation Liquid. In 100mM TRIS-HCl, pH 7.5, containing 14mM β -mercaptoethanol, 0.5mM EDTA, 0.5mM PMSF and 10% glycerol.

Purity $\geq 99\%$ (SDS-PAGE)

Purity Detail Affinity purified.

Source Produced in Sf9 insect cells. Produced in a baculovirus expression system.

Specific Activity $\geq 600\text{U/mg}$ protein. One unit is defined as the amount of enzyme that synthesizes 1nmol of poly(ADP-ribose) per min. at 25°C, pH 7.5.

UniProt ID P09874

Last modified: May 29, 2024



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)