

PARP-2 (mouse), (recombinant) (high purity)

Citations: 5

[View Online »](#)

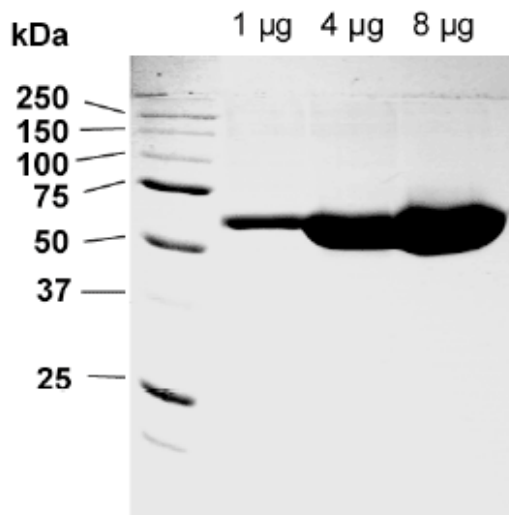
Ordering Information

[Order Online »](#)

ALX-201-0640	20µg
--------------	------

Manuals, SDS & CofA

[View Online »](#)



Quality Control: 12% 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 2
Concentration	~1mg/ml
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 1mM DTT, 0.5mM EDTA, 0.5mM PMSF and 20% glycerol.
Purity	≥98% (SDS-PAGE)
Purity Detail	Affinity purified.
Source	Produced in Sf9 insect cells (full length). Produced in a baculovirus expression system.
Specific Activity	≥200U/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. See also separate protocol (link below).
UniProt ID	O88554

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com