

PARP-3 (human), (recombinant) (high purity) (His-tag)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-252-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Poly(ADP-ribose) polymerase 3
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 100mM sodium chloride, 50mM imidazole, 0.2% NP-40 and 10% glycerol.
MW	~60kDa
Purity	≥98% (SDS-PAGE)
Source	Produced in Sf21 insect cells. Human full-length PARP-3 is fused to a C-terminal His-tag. Produced in a baculovirus expression system.
UniProt ID	Q9Y6F1

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)