

# PARP-1 (human), (recombinant) (high purity) (His-tag)

Citations: 3

[View Online »](#)

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 1
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 100mM sodium chloride, 50mM imidazole, 0.2% NP-40 and 10% glycerol.
Purity	≥98% (SDS-PAGE)
Source	Produced in Sf21 insect cells. Human full-length PARP-1 is fused to a Myc-tag and a His-tag. Produced in a baculovirus expression system.
Specific Activity	≥1'200U/mg protein.
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)