

PARP-1 (catalytic domain) (human), (recombinant) (His-tag)

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 10mM TRIS-HCl, pH 7.5, containing 300mM sodium chloride, 250mM imidazole, 2mM β -mercaptoethanol, 0.2% NP-40 and 10% glycerol.
MW	~46kDa
Purity	≥98% (SDS-PAGE)
Source	Produced in Sf21 insect cells. Human PARP-1 (656-1014) with Myc-tag and HIS-tag. Produced in a baculovirus expression system.
UniProt ID	P09874

Last modified: June 11, 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