

PRODUCT INFORMATION



PARP1 (human, recombinant)

Item No. 32561

Overview and Properties

Synonyms: ADP-ribose Transferase 1, ADPRT, PARP1, PARP-1, NAD⁺ADP-ribosyltransferase 1, Poly(ADP-ribose) Polymerase 1

Source: Active recombinant human N-terminal GST-tagged PARP1 expressed in insect cells

Amino Acids: 2-1,041 (full length)

Uniprot No.: P09874

Molecular Weight: 139 kDa

Storage: -80°C (as supplied)

Stability: ≥6 months

Purity: **batch specific** (≥80% estimated by SDS-PAGE)

Supplied in: 40 mM Tris, pH 8.0, with 110 mM sodium chloride, 2.2 mM potassium chloride, 0.5 mM glutathione, 3 mM DTT, and 20% glycerol

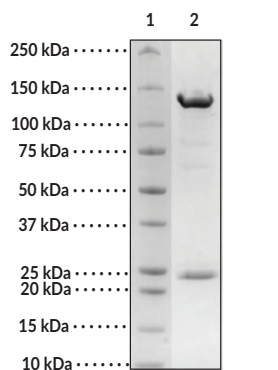
Protein

Concentration: **batch specific** mg/ml

Activity: Ribosylation activity measured by chemiluminescent assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

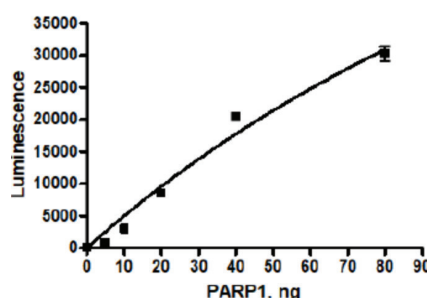
Images



Lane 1: MW Markers
Lane 2: PARP1 (2 µg)

SDS-PAGE Analysis of PARP1.

Representative gel image shown; actual purity may vary between each batch.



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/11/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Poly(ADP-ribose) polymerase 1 (PARP1) is an ADP-ribosylating enzyme that has roles in DNA repair, maintenance of genomic integrity, and transcriptional regulation.^{1,2} It is composed of an N-terminal DNA-binding domain (DBD) that contains the nuclear localization signal (NLS) and three zinc fingers that mediate PARP1 self-assembly and activation, a central automodification domain, and a highly conserved C-terminal catalytic domain.^{1,3} PARP1 is ubiquitously expressed and localizes to the nucleus where it is recruited to sites of DNA damage induced by a variety of cellular stressors, including genomic, oxidative, inflammatory, or metabolic stress.^{1,4} Binding of PARP1 to DNA single- or double-strand breaks activates its poly(ADP-ribosylation (PARylation) catalytic activity. PARP1 PARylates itself, increasing its activity and recruiting additional DNA repair proteins to sites of damaged DNA, as well as PARylates other proteins, including transcription factors, to facilitate DNA damage repair.^{1,4} PARP1 is subject to additional PTMs, including phosphorylation, methylation, and acetylation, that regulate its catalytic and DNA-binding activities.⁵ Increased PARP1 levels have been found in tumors isolated from patients with a variety of cancers, including soft tissue sarcoma and ovarian or squamous cell carcinomas, and are associated with decreased survival. Cayman's PARP1 (human, recombinant) protein can be used for enzyme assay applications. This protein consists of 1,040 amino acids and has a calculated molecular weight of 139 kDa.

References

1. Chaudhuri, A.R. and Nussenzweig, A. The multifaceted roles of PARP1 in DNA repair and chromatin remodelling. *Nat. Rev. Mol. Cell Biol.* **18**(10), 610-621 (2017).
2. Ko, H.L. and Ren, E.C. Functional aspects of PARP1 in DNA repair and transcription. *Biomolecules* **2**(4), 524-548 (2012).
3. Alemasova, E.E. and Lavrik, O.I. Poly(ADP-ribosyl)ation by PARP1: Reaction mechanism and regulatory proteins. *Nucleic Acids Res.* **47**(8), 3811-3827 (2019).
4. Luo, X. and Kraus, W.L. On PAR with PARP: Cellular stress signaling through poly(ADP-ribose) and PARP-1. *Genes Dev.* **26**(5), 417-432 (2012).
5. Pazzaglia, S. and Pioli, C. Multifaceted role of PARP-1 in DNA repair and inflammation: Pathological and therapeutic implications in cancer and non-cancer diseases. *Cells* **9**(1), 41 (2020).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM