

Recombinant PARP-1 protein

Catalog No: 31238

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.25 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.25 µg/µl in 20 mM Tris-HCl, pH 7.3, 125 mM NaCl, 0.2 mM EDTA and 20% glycerol.

Background: PARP-1, poly (ADP-ribose) polymerase family, member 1, (also known as PARP1, ADPRT, ADP-ribosyltransferase NAD⁺, or PPOL, poly(ADP-ribose) polymerase) is an abundant and ubiquitous nuclear enzyme involved in cellular processes such as DNA damage repair, the regulation of chromatin structure and the control of gene expression. PARP-1 catalyzes the poly(ADP-ribosyl)ation of a variety of acceptor proteins involved in chromatin architecture and in DNA metabolism, including core histones, the linker histone H1 and a variety of transcription factors.

For additional information on PARP-1, please see the review article PARP-1: An Abundant and Ubiquitous Protein with Roles in Many Cellular Processes in the Targets & Applications section or our website.

Protein Details: Recombinant PARP-1 protein was expressed in Sf9 insect cells as the full-length human protein (accession number NP_001609) with an N-terminal FLAG tag. The molecular weight of the protein is ~115 kDa.

Application Notes: Recombinant PARP-1 protein is suitable for use in assays that modify histones and CTCF by poly ADP-ribosylation. A recommended starting point is 100 ng recombinant protein per reaction.

Storage and Guarantee: This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.