
PARP1 Protein

(Catalog No. E53000)

Description

Human PARP1 recombinant protein, 662-1014aa

Molecular Weight (kDa)

39.6 kDa

Formulation

20 mM Tris buffer(pH 8.0) containing 10% glycerol 1mM DTT.

Source

Expressed in E.coli expression system

Size

25 µg

Storage Temp & Conditions

Stored at 4°C short term (1-2 weeks). Store at -20°C or -80°C long term

Purification Notes

Purity :> 95% by SDS-PAGE

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array

Ordering Information:

Products	Size	Cat. No.
PARP1 Protein	25 µg	E53000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.