
PARP2 Protein

(Catalog No. E54000)

Description

Human PARP2 (GenBank Accession No.NM_005484), amino-acids 2-583 with N-terminalGST tag

Molecular Weight (kDa)

92 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 Mm KCl, 18 mMglutathione, 10% glycerol, and 3 mM DTT.

Source

Produced in Sf21 cells.

Size

10 µg

Storage Temp & Conditions

-80°C. After opening, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.

Purification Notes

Purity: ≥80% (SDS-PAGE)

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
PARP2 Protein	10 µg	E54000-1

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