
PARP3 Protein

(Catalog No. E54001)

Description

Human PARP3, long isoform (GenBank Accession No. NM_001003931), full length with N- terminal GST tag.

Molecular Weight (kDa)

87 kDa

Formulation

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

Source

Expressed in a Baculovirus infected Sf9 cell expression system.

Size

10 µg

Storage Temp & Conditions

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

Purification Notes

Purity: ≥50% (SDS-PAGE)

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
PARP3 Protein	10 µg	E54001-1

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