
XRCC3 Protein

(Catalog No. E72012)

Description

Human XRCC3 full-length ORF (AAH01036.1, 1 a.a. - 346 a.a.) recombinant protein with GST-tag at N-terminal.

Molecular Weight (kDa)

63.69 kDa

Formulation

50 mM Tris-HCl, 10 mM reduced Glutathione, pH=8.0 in the elution buffer.

Source

In vitro wheat germ expression system

Size

10 µg

Storage Temp & Conditions

Store at -80°C. Aliquot to avoid repeated freezing and thawing. Best use within three months from the date of receipt of this protein.

Purification Notes

Glutathione Sepharose 4 Fast Flow

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
XRCC3 Protein	10 µg	E72012-1

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