

---

## NF-kb p50 Protein

(Catalog No. E62001)

---

### Description

Human NFKB1 full-length ORF ( AAH51765.1, 1 a.a. - 969 a.a.) recombinant protein with GST-tag at N-terminal.

### Molecular Weight (kDa)

131.8 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.

### 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.
NF-kb p50 Protein	10 µg	E62001 -1

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