

---

## MIZF Protein

(Catalog No. E42005)

---

### Description

Human MIZF partial ORF ( NP\_056332, 1 a.a. - 110 a.a.) recombinant protein with GST-tag at N-terminal

### Molecular Weight (kDa)

37.84 kDa

### Formulation

51 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.
MIZF Protein	10 µg	E42005-1

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