
c-Myc Protein

(Catalog No. E62000)

Description

Human MYC full-length ORF (ENSP00000259523, 1 a.a. - 439 a.a.) recombinant protein with GST-tag at N-terminal

Molecular Weight (kDa)

75.2 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.
c-Myc Protein	10 µg	E62000 -1

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