## c-Myc (human) Recombinant monoclonal antibody (9E10) (Mouse F(ab)2)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS458-0100

100µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Use/Stability** Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Proto-oncogene c-Myc, Transcription factor p64, Class E

basic helix-loop-helix protein 39, BHLHE39

**Application** IF, IHC (FS), IHC (PS), IP, WB

**Clone** 9E10

**Formulation** Liquid. In PBS containing 0.02% Proclin 300.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to aa 408-439 from C-

terminus of human c-myc.

Isotype F(ab)2

Purity Detail Immobilized metal affinity chromatography purified.

Species Reactivity Human

**Specificity** Recognizes epitope sequence EQKLISEEDL, located in

aa 410-419 of human c-myc protein.

**Technical Info / Product Notes**This is a recombinant mouse Fab2 fragment, based on the

original mouse IgG1 format, created for improved compatibility with existing reagents, assays and

techniques.

P01106

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



eu@enzolifesciences.com