

# c-Myc (human) Recombinant monoclonal antibody (9E10) (Mouse IgG2bk)

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                           |   |
|---------------------------|---|
| <b>Use/Stability</b>      | Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C. |
| <b>Handling</b>           | Avoid freeze/thaw cycles.   |
| <b>Short Term Storage</b> | +4°C  |
| <b>Long Term Storage</b>  | -20°C   |
| <b>Shipping</b>           | Blue Ice  |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                                       |  |
|---------------------------------------|--|
| <b>Alternative Name</b>               | Proto-oncogene c-Myc, Transcription factor p64, Class E basic helix-loop-helix protein 39, BHLHE39   |
| <b>Application</b>                    | IF, IHC (FS), IHC (PS), IP, WB   |
| <b>Clone</b>                          | 9E10   |
| <b>Formulation</b>                    | Liquid. In PBS containing 0.02% Proclin 300.   |
| <b>Host</b>                           | Mouse  |
| <b>Immunogen</b>                      | Synthetic peptide corresponding to aa 408-439 from C-terminus of human c-myc.  |
| <b>Isotype</b>                        | IgG2bk   |
| <b>Purity Detail</b>                  | Protein A affinity purified.   |
| <b>Species Reactivity</b>             | Human  |
| <b>Specificity</b>                    | Recognizes epitope sequence EQKLISEEDL, located in aa 410-419 of human c-myc protein.  |
| <b>Technical Info / Product Notes</b> | This is a reformatted mouse IgG2b antibody, based on the original mouse IgG1 format, created for improved compatibility with existing reagents, assays and techniques. |

UniProt ID

P01106

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)