

# c-Myc (human) Recombinant monoclonal antibody (9E10) (Mouse Fab fragment $\kappa$ ) (His-tag)

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

Ordering Information

[Order Online »](#)

ENZ-ABS459-0100	100 $\mu$ g
-----------------	-------------

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>	Fab fragment κ
<b>Purity Detail</b>	Immobilized metal affinity chromatography 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 recombinant mouse Fab fragment, based on the original mouse IgG1 format, created for improved compatibility with existing reagents, assays and techniques.
<b>UniProt ID</b>	P01106

Worry-free Guarantee

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



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)