

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

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

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	Rat
Immunogen	Synthetic peptide corresponding to aa 408-439 from C-terminus of human c-myc.
Isotype	IgG2bk
Purity Detail	Protein A affinity 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 chimeric rat 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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