c-Myc (human) Recombinant monoclonal antibody (9E10) (Human IgG4κ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS474-0200

200µg

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 Human

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

terminus of human c-myc.

lsotype lgG4k

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 NotesThis is a chimeric human antibody, 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