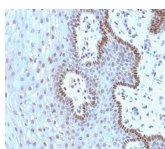


c-Myc (EQKLISEEDL) Monoclonal Mouse Antibody (9E10.3)



Product Description

The epitope of this antibody spans the Myc tag epitope, aa 410-419 (EQKLISEEDL), which is a specific portion of an alpha helical region of human c-myc protein. Because the epitope is located at the dimerization site of the protein, the antibody is not sensitive for detecting endogenous c-myc, making it suitable for detecting Myc tagged proteins or myc overexpression. This MAb shows no cross-reaction with v-myc. c-myc is involved in the control of cell proliferation and differentiation and is amplified and/or overexpressed in a variety of tumors. Over-expression of c-myc protein occurs frequently in luminal cells of prostate intraepithelial neoplasia as well as in most primary carcinomas and metastatic disease.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0596, Anti-c-Myc (9E10.3)

Product attributes

Antibody number	#0596
Antibody reactivity (target)	c-Myc
Antibody type	Primary, Tag Antibody
Host species	Mouse
Clonality	Monoclonal
Clone	9E10.3
Isotype	IgG1, kappa
Molecular weight	62-64 kDa
Synonyms	Class E basic helix-loop-helix protein 39 (bHLHe39); MRTL; Myc2; Nlard; Nird; Proto-oncogene c-Myc; RNCMYC; Transcription factor p64; Transcriptional regulator Myc-A; V-Myc avian myelocytomatosis viral oncogene homolog
Human gene symbol	MYC
Entrez gene ID	4609
SwissProt	P01106
Unigene	202453
Immunogen	A synthetic peptide, corresponding to aa 408-439 (AEEQKLISEEDLLRKRREQLKHKLEQLRNSCA) from C-terminus of human c-myc, coupled to KLH.
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	Immunohistochemistry formalin-fixed 1-2 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Immunofluorescence 1-2 ug/mL. Does not react with mouse, rat, or chicken, others not tested. Optimal dilution for a specific application should be determined by user
Positive control	HL-60 cells or breast carcinoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Cancer, Cell cycle, Transcription factors
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 21, 2024 at 07:55:02 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-c-myc-9e10/>