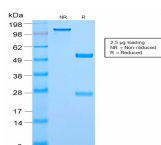


# c-Myc (EQKLISEEDL) Recombinant Monoclonal Rabbit Antibody (MYC2895R)



## 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2895, Anti-c-Myc Oncoprotein (MYC2895R)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

<b>Antibody number</b>	#2895
<b>Antibody reactivity (target)</b>	c-Myc Oncoprotein
<b>Antibody type</b>	Primary
<b>Host species</b>	Rabbit
<b>Clonality</b>	Monoclonal, Recombinant Monoclonal
<b>Clone</b>	MYC2895R
<b>Isotype</b>	IgG
<b>Molecular weight of antigen</b>	62-64 kDa
<b>Synonyms</b>	Class E basic helix-loop-helix protein 39 (bHLHe39); MRTL; Myc2; Niard; Nird; Proto-oncogene c-Myc; RNCMYC; Transcription factor p64; Transcriptional regulator Myc-A; V-Myc avian myelocytomatosis viral oncogene homolog
<b>Human gene symbol</b>	MYC
<b>Entrez gene ID</b>	4609
<b>SwissProt</b>	P01106
<b>Unigene</b>	202453
<b>Immunogen</b>	A synthetic peptide, corresponding to aa 408-439 (AEEQKLISEEDLLRKRREQLKHKLEQLRNSCA) from C-terminus of human c-myc, coupled to KLH.
<b>Verified antibody applications</b>	Flow (intracellular) (verified)
<b>Antibody target cellular localization</b>	Nucleus
<b>Species reactivity</b>	Human
<b>Positive control</b>	K562, HL-60 cells. Cervical Carcinoma.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Antibody research areas</b>	Cancer, Cell cycle, Transcription factors

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 June 3, 2026 at 03:41:25 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/c-myc-recombinant-monoclonal-rabbit-antibody-myc2895r/>