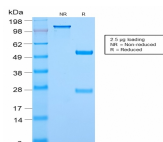


# c-Myc (EQKLISEEDL) Recombinant Monoclonal Rabbit Antibody (MYC2895R)



## 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) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

## Product attributes

Antibody number	#2895
Antibody reactivity (target)	c-Myc Oncoprotein
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	MYC2895R
Isotype	IgG
Molecular weight	62-64 kDa
Synonyms	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
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	Flow (intracellular) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Positive control	K562, HL-60 cells. Cervical Carcinoma.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	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.
Antibody research areas	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 August 31, 2025 at 02:04:55 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/c-myc-recombinant-monoclonal-rabbit-antibody-myc2895r/>