## Biotin Recombinant Monoclonal Rabbit Antibody (BTN/2032R)



## **Product Description**

This antibody recognizes both the free and protein-conjugated (either soluble or cell bound) form of biotin. This MAb is highly specific to biotin and shows no cross-reaction with other structurally related compounds. It has a very high affinity for biotin and is excellent for use in various amplification techniques. In some applications, localization of biotinylated probes with avidin produces unacceptably high background staining. Anti-biotin antibody may be substituted to decrease this noise.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2032, Anti-Biotin (BTN/2032R)

## Product attributes

Product attributes		
Antibody number	#2032	
Antibody reactivity (target)	Biotin	
Antibody type	Primary, Tag Antibody	
Host species	Rabbit	
Clonality	Recombinant Monoclonal	
Clone	BTN/2032R	
Isotype	IgG	
Molecular weight	244Da	
Synonyms	Vitamin B7 or Vitamin H	
Human gene symbol	Not Applicable	
Entrez gene ID	Not Applicable	
SwissProt	ot Applicable	
Unigene	Not Applicable	
Immunogen	Biotinylated sheep immunoglobulin	
Verified antibody applications	HC (FFPE) (verified)	
Species reactivity	All species	
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Immunohistology (formalin): 0.5-1 ug/mL for 30 minutes at RT, In situ Nucleic Acid Hybridization, No special pretreatment is required for the immunohistochemical staining of formalin-fixed paraffirin-embedded tissues., Western blotting, Optimal dilution for a specific application should be determined by user	
Positive control	Biotinylated proteins in solution or on tissues	
Positive control Shipping condition	, ,,	
	Biotinylated proteins in solution or on tissues	
Shipping condition	Biotinylated proteins in solution or on tissues Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,	
Shipping condition Storage Conditions	Biotinylated proteins in solution or on tissues  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when	
Shipping condition Storage Conditions Shelf life	Biotinylated proteins in solution or on tissues  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended	

Call us: 800-304-5357 Email: btinfo@biotium.com

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,	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 8, 2025 at 04:07:12 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-recombinant-monoclonal-rabbit-antibody-btn-2032r/">https://biotium.com/product/biotin-recombinant-monoclonal-rabbit-antibody-btn-2032r/</a>