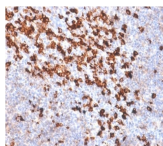


Biotin Recombinant Monoclonal Rabbit Antibody (BTN/2032R)



Product Description

This antibody recognizes both the free and protein-conjugated (either soluble or cell bound) form of biotin. This MAbs 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 order@biotium.com 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

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	Not Applicable
Unigene	Not Applicable
Immunogen	Biotinylated sheep immunoglobulin
Verified antibody applications	IHC (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, Immunohistochemistry (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 paraffin-embedded tissues., Western blotting, Optimal dilution for a specific application should be determined by user
Positive control	Biotinylated proteins in solution or on tissues
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
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 # 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 September 17, 2025 at 04:13:53 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-recombinant-monoclonal-rabbit-antibody-btn-2032r/>