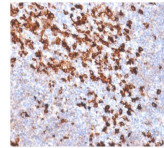


## Biotin Recombinant Monoclonal Rabbit Antibody (BTN/2032R)



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

### 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 [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 2032, Anti-Biotin (BTN/2032R)**

Antibody #	prefix	Conjugation	Ex/Em	Concentration	Storage Buffer
BNC04	CF@405S		404/431 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC05	CF@405M		408/452 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC88	CF@488A		490/515 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC43	CF@543		541/560 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC55	CF@555		555/565 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC68	CF@568		562/583 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC94	CF@594		593/614 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC40	CF@640R		642/662 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC47	CF@647		650/665 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC61	CF@660R		663/682 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC80	CF@680		681/698 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC81	CF@680R		680/701 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC70	CF@770		770/797 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCR	R-PE (PE)		496, 546, 565/578 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCA	APC		650/660 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCP	PerCP		482/677 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCB	Biotin	N/A	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCAP	Alkaline Phosphatase	N/A	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCH	Horseradish Peroxidase	N/A	N/A	0.1 mg/mL	PBS, 0.05% BSA, no azide
BNUB	Purified, with BSA	N/A	N/A	0.2 mg/mL	PBS, 0.05% BSA, 0.05% azide
BNUM	Purified, BSA-free	N/A	N/A	1 mg/mL	PBS, no BSA, no azide

### References

Uccellini MB, et al. 2008. J. Immunol. 181:5875.[Green NM, et al. 2012. J Biol Chem. 287:3978.

### Product attributes

Antibody number	2032
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
Species reactivity	All species
Applications	Immunofluorescence, Immunohistology (formalin), Flow cytometry, Western, In situ hybridization
Application notes	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 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)

This datasheet was generated on September 28, 2020 at 11:29:01 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-recombinant-monoclonal-rabbit-antibody-btn-2032r/>