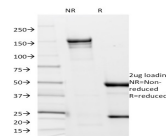


Borrelia burgdorferi p41 Flagellin Monoclonal Mouse Antibody (6802)



Product Description

This antibody recognizes the p41 flagellar antigen (Flagellin) of the bacteria *Borrelia burgdorferi*. *Borrelia burgdorferi* is a species of gram-negative bacteria of the spirochete class of the genus *Borrelia*. *B. burgdorferi* resembles other spirochetes in that it is a highly specialized, motile, two-membrane, spiral shaped bacteria which lives primarily as an extracellular pathogen. *Borrelia burgdorferi* is a primary cause of Lyme disease, a vector-borne, multisystem inflammatory disease which is transmitted to humans by the bite of ticks of the *Ixodes ricinus* complex.

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 0144, *Borrelia burgdorferi* p41 Flagellin Monoclonal Mouse Antibody (6802)

Product attributes

Antibody number	#0144
Antibody reactivity (target)	<i>Borrelia burgdorferi</i> p41 Flagellin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	6802
Isotype	IgG2a, kappa
Molecular weight	~41 kDa
Synonyms	<i>B. burgdorferi</i> ; Flagellar filament 41kDa core protein; Lyme disease
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	A crude sonicate of <i>Borrelia</i> spirochetes
Species reactivity	<i>Borrelia burgdorferi</i>
Antibody application notes	For coating for ELISA, order Ab without BSA. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	<i>Borrelia burgdorferi</i> extract or infected cells. Tissue.
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
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
Antibody research areas	Microbiology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), 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 January 1, 2026 at 01:26:06 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/borrelia-burgdorferi-p41-flagellin-monoclonal-mouse-antibody-6802/>