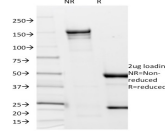


Borrelia burgdorferi p41 Flagellin Monoclonal Mouse Antibody (6802)



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

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
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, LI-COR Bioscience.

This datasheet was generated on July 17, 2024 at 09:33:32 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/borrelia-burgdorferi-p41-flagellin-monoclonal-mouse-antibody-6802/>