Pseudomonas Aeruginosa Serotype 5C Monoclonal Mouse Antibody (320/65)

Product Description

This antibody is specific for serotype 5C and does not react with other species. Pseudomonas aeruginosa is a Gram-negative, aerobic, rod-shaped bacterium with unipolar motility. It is an opportunistic pathogen of plants and humans and can infect the urinary tract, respiratory and gastrointestinal system, soft tissues, bones and joints leading to severe systemic infections of immunosuppressed patients in hospitals. P. aeruginosa secretes a variety of pigments, including pyocyanin (blue-green), fluorescein (pyoverdin), and pyorubin (red-brown). This organism can achieve anaerobic growth with nitrate as a terminal electron acceptor, and, in its absence, it is also able to ferment arginine by substrate-level phosphorylation. Adaptation to microaerobic or anaerobic environments is essential for certain lifestyles of P. aeruginosa, such as during lung infection in cystic fibrosis patients where thick layers of alginate surrounding bacterial mucoid cells can limit the diffusion of oxygen.

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 0239, Anti-Pseudomonas Aeruginosa Serotype 5C (320/65)

Product attributes

Call us: 800-304-5357

i ioddot attiibates			
Antibody number	#0239		
Antibody reactivity (target)	Pseudomonas Aeruginosa Serotype 5C		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	320/65		
Isotype	IgM, kappa		
Molecular weight	Not Known		
Synonyms	Pseudomonas aeruginosa serotype 5C		
Human gene symbol	Not Applicable		
Entrez gene ID	Not Applicable		
SwissProt	Not Applicable		
Unigene	Not Applicable		
Immunogen	Pseudomonas aeruginosa serotype 5C		
Species reactivity	Pseudomonas aeruginosa serotype 5C		
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	Pseudomonas aeruginosa serotype 5C 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		
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.		

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, 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 November 2, 2025 at 03:58:58 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/pseudomonas-aeruginosa-serotype-5c-monoclonal-mouse-antibody-320-65/