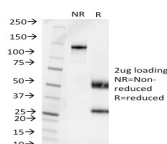


Pseudomonas Aeruginosa Serotype 6C Monoclonal Mouse Antibody (1200/472)



Product Description

This antibody is specific for serotype 6C and does not react with other species. *Pseudomonas aeruginosa* is Gram-negative, aerobic, rod-shaped bacteria with unipolar motility. An opportunistic human pathogen, *P. aeruginosa* is also an opportunistic pathogen of plants. *P. aeruginosa* bacteria are clinically important because they are resistant to most antibiotics and they are capable of surviving in conditions that few other organisms can tolerate. *Pseudomonas* is often encountered in hospital and clinical work because it is a major cause of hospital acquired (nosocomial) infections. Its main targets are immuno-compromised individuals, burn victims, and individuals on respirators or with indwelling catheters. Additionally, these pathogens colonize the lungs of cystic fibrosis patients. *P. aeruginosa* is often identified by its pearlescent appearance and grape-like odor in vitro. Definitive clinical identification of *P. aeruginosa* includes identifying the production of both pyocyanin and fluorescein as well as its ability to grow at 42C. *P. aeruginosa* is capable of growth in diesel and jet fuel, where it is known as hydrocarbon utilizing microorganisms (or HUM bugs), causing microbial corrosion.

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 0240, Anti-Pseudomonas Aeruginosa Serotype 6C (1200/472)

Product attributes

Antibody number	#0240
Antibody reactivity (target)	Pseudomonas Aeruginosa Serotype 6C
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	1200/472
Isotype	IgG1, kappa
Molecular weight	Not Known
Synonyms	Pseudomonas aeruginosa serotype 6C
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Pseudomonas aeruginosa serotype 6C
Species reactivity	Pseudomonas aeruginosa serotype 6C
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 6C 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 (<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 December 12, 2025 at 06:36:56 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/pseudomonas-aeruginosa-serotype-6c-monoclonal-mouse-antibody-1200-472/>