



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Campylobacter jejuni Monoclonal Mouse Antibody (2E-10)

Product Description

This MAb is specific for a non-flagellar antigen of both C. jejuni Types 1 and 2. It reacts with a soluble excreted antigen in EIA. This determinant is unaffected by frozen storage of specimens, unlike antibodies to flagellar antigens which require fresh cultured organisms. Campylobacter jejuni is a Gram-negative, microaerophilic, slender, flagellate, spiral bacterium. It is the major reported cause of bacterial foodborne infection in the United States and is also associated with Guillian-Barre syndrome. Campylobacteriosis is an infectious disease caused by bacteria of the genus Campylobacter. The Campylobacter organism is actually a group of spiral shaped bacteria that can cause disease in humans and animals. Most human illness is caused by one species, called Campylobacter jejuni, but 1% of human Campylobacter cases are caused by other species.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Catalog number key for antibody number 0434, Anti-Campylobacter jejuni (2E-10.)

Product attributes

Antibody number	#0434
Antibody reactivity (target)	Campylobacter jejuni
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	2E-10.
Isotype	IgM, kappa
Molecular weight	Not Known
Synonyms	Campylobacter jejuni
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Total sonicate of Campylobacter jejuni
Species reactivity	Campylobacter jejuni
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	Campylobacter jejuni infected stomach biopsy
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, 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)
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
Antibody research areas	Infectious disease, Microbiology
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.

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 14, 2026 at 12:42:28 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mouse-monoclonal-anti-campylobacter-jejuni-2e-10/>