



Glowing products for science

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

E. coli O157 Monoclonal Mouse Antibody (9.88)

Product Description

This MAb shows specificity to E.coli O157 in a simple ELISA. Escherichia coli are Gram negative bacterium that are commonly found in the lower intestine of warm-blooded organisms (endotherms). Their serological types are determined in combination with somatic antigens (O group: O1-O173) and flagella antigens (H type: H1-H56). The E. coli that cause intestinal infectious diseases including diarrhea, acute gastritis or colitis are referred to as pathogenic E. coli, which are classified into the following 4 groups according to differences in the mode of pathogenicity; enteropathogenic E. coli (EPEC), enteroinvasive E. coli (EIEC), enterotoxigenic E. coli (ETEC) and enterohemorrhagic E. coli (EHEC). Although the identification of pathogenic E. coli requires verification of their pathogenicity, pathogenic E. coli often have specific serotypes; therefore, typing of the serogroup and serotype is necessary in screening pathogenic E. coli.

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 0207, Anti-E. coli O157 (9.88)**

Product attributes

Antibody number	#0207
Antibody reactivity (target)	E. coli O157
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	9.88
Isotype	IgM, kappa
Molecular weight	Not Known
Synonyms	E. coli O157
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	A crude sonicate of E. coli O157
Species reactivity	E. coli O157
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	E. coli O157 extract or infected cells or tissue
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 31, 2026 at 08:09:11 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mouse-monoclonal-anti-e-coli-o157-9-88/>