

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.

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

This datasheet was generated on July 17, 2024 at 09:20:42 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mouse-monoclonal-anti-e-coli-o157-9-88/>