

Campylobacter jejuni Monoclonal Mouse Antibody (2E-10.)

This MAb is specific for a non-flagellar antigen of both C. jejuni Types 1 and 2.

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.



Product attributes

Antibody number	0434
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
Applications	Immunofluorescence, ELISA
Application notes	ELISA, for coating order Ab without BSA and use at 1-5 ug/mL, Immunofluorescence 1-2 ug/mL, Optimal dilution for a specific application should be determined by user
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)
Research areas	Infectious disease, Microbiology