

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.

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 <i>Campylobacter jejuni</i>
Species reactivity	<i>Campylobacter jejuni</i>
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	<i>Campylobacter jejuni</i> 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