

Lewis-Y Monoclonal Mouse Antibody (A70-C/C8)

Product Description

This antibody recognizes a carbohydrate epitope common to the tumor-associated Lewis Y and Lewis b antigens (Fucalpha1-2Galbeta1-4/3[Fucalpha1-3/4]GlcNAcbeta). Its specificity was established without doubt with a panel of 86 synthetic mono- and oligosaccharidic structures. Lewis Y is expressed in large bowel tumors and colorectal carcinomas. It may be useful in the classification of human renal and bladder tumors. The Lewis Y antigen has been evaluated as a clinical marker for the diagnosis and prognosis of cholangiocarcinoma, hepatocellular carcinoma and breast cancer.

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

Product attributes

Antibody number	#1411
Antibody reactivity (target)	Lewis-Y
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	A70-C/C8
Isotype	IgM
Molecular weight	Multiple
Synonyms	Lewis Y antigen
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
UniGene	Not Applicable
Immunogen	Live Ls174T cells (human colon carcinoma cell line)
Antibody target cellular localization	Cytoplasmic, Plasma membrane, Secreted (extracellular)
Species reactivity	Human
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	Lung, Colon, Bladder or Hepatocellular Carcinoma
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	Cancer
Tumor expression	Bladder cancer, Colorectal cancer, Renal cancer

This datasheet was generated on August 29, 2024 at 08:29:39 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-lewis-y-a70-cc8/>