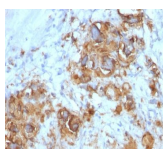


CA19-9 / Sialyl-Lewis A Monoclonal Mouse Antibody (121SLE)



Product Description

CA19-9, a carbohydrate epitope expressed on a high MW (>400 kDa) mucin glycoprotein, is a sialyl Lewisia structure which is synthesized from type 1 blood group precursor chains. It is present in individuals expressing the Lewisia and/or Lewisb blood group antigens. In normal tissues, sialyl Lewisia antigen is present in ductal epithelium of the breast, kidney, salivary gland, and sweat glands. Its expression is greatly enhanced in serum as well as in the majority of tumor cells in gastrointestinal (GI) carcinomas, including adenocarcinomas of the stomach, intestine, and pancreas. Preoperative elevated CA19-9 levels in patients with stage I pancreatic carcinoma decrease to normal values following surgery. When used serially, CA19-9 can predict recurrence of disease prior to radiographic or clinical findings. This MAb is excellent for staining of formalin-fixed, paraffin-embedded tissues.

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 0025, Anti-CA19-9|Sialyl Lewis A (121SLE)

Product attributes

Antibody number	#0025
Antibody reactivity (target)	CA19-9, Sialyl Lewis A
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	121SLE
Isotype	IgM, kappa
Molecular weight	>400 kDa
Synonyms	CA19.9, Ovarian tumor antigen, Pancreatic tumor antigen, Sialyl Lewis a
Human gene symbol	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Precipitin lines obtained after immuno-diffusion using MAb 116-NS-19-9 and mucins isolated from an ovarian cyst of a Lewis A B- patient (0Le).
Antibody target cellular localization	Secreted (extracellular), Vesicular
Cell/tissue expression	Epithelial cells
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Antibody application notes	Immunohistology (formalin), No special pretreatment is required for the immunohistochemical staining of formalin-fixed, paraffin-embedded tissues, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL
Positive control	Stomach or Colon 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, Hematology
Tumor expression	Gastrointestinal cancer, Pancreatic cancer
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 September 13, 2025 at 07:55:18 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-ca19-9-sialyl-lewis-a-121sle/>