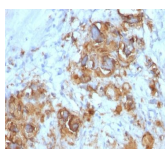


# CA19-9 / Sialyl-Lewis A Monoclonal Mouse Antibody (121SLE)



## 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

<b>Antibody number</b>	#0025
<b>Antibody reactivity (target)</b>	CA19-9, Sialyl Lewis A
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	121SLE
<b>Isotype</b>	IgM, kappa
<b>Molecular weight</b>	>400 kDa
<b>Synonyms</b>	CA19.9, Ovarian tumor antigen, Pancreatic tumor antigen, Sialyl Lewis a
<b>Human gene symbol</b>	Not Known
<b>Entrez gene ID</b>	Not Known
<b>SwissProt</b>	Not Known
<b>Unigene</b>	Not Known
<b>Immunogen</b>	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).
<b>Antibody target cellular localization</b>	Secreted (extracellular), Vesicular
<b>Cell/tissue expression</b>	Epithelial cells
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	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
<b>Positive control</b>	Stomach or Colon Carcinoma
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer, Hematology
<b>Tumor expression</b>	Gastrointestinal cancer, Pancreatic cancer
<b>Product origin</b>	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	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
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 November 23, 2024 at 10:59:50 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/>