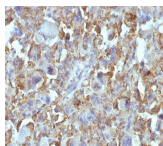


Small Cell Lung Cancer Monoclonal Mouse Antibody (MOC-52)



Product Description

This MAb reacts with a membrane-associated protein present in normal and malignant neuroendocrine tissues including small cell lung cancer (SCLC). It stains neural and a variable number of endocrine tissues and in the lung it reacts preferentially with SCLC and carcinoids. Its epitope is destroyed during formalin fixation. This antibody was categorized during the First International Workshop on Small Cell Lung Cancer Antigens held in London in April 1987. There are two major types of Lung Carcinoma: non-small cell, which accounts for 80% of all cases; and small cell, which accounts for roughly 20% of all lung cancers reported. The lung continues to be a customary place for cancer migration from tumors elsewhere in the body. Treatment depends on the specific cell type of the cancer, level of progression and status of the individual patient.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0329, Anti-Small Cell Lung Cancer (MOC-52)

Product attributes

Antibody number	#0329
Antibody reactivity (target)	Small Cell Lung Cancer
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MOC-52
Isotype	IgG1, kappa
Molecular weight	145 kDa
Synonyms	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Lung Cancer-associated antigen isolated from small cell lung carcinoma-derived cell line
Antibody target cellular localization	Plasma membrane
Expected antibody applications	IHC (frozen) (published for clone)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL, Immunocytochemistry frozen 0.5-1.0 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	Human small cell lung carcinoma, adrenal gland
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, 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
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.
Tumor expression	Lung cancer

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Souhami RL, Beverly PCL, Bobrow LG. (1987) Antigens of small-cell lung cancer: first international workshop. Lancet 2:325-326. (clone characterization)
2. J Clin Oncol (1989) 7(11): 1614-1620. (IHC, frozen)

This datasheet was generated on July 12, 2025 at 02:47:06 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-small-cell-lung-cancer-sclc-moc-52/>