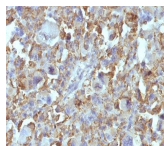


Histiocytoma Marker Monoclonal Mouse Antibody (D11)



Product Description

In Western blotting, this antibody detects an antigen of 125 kDa in human liver and 135 kDa in tumors of histiocytic origin. Comparative study of this MAb and a standard CD68 MAb showed that their antigens are different. Its antigen in all macrophage types studied is located on the plasma membrane and within cytoplasmic structures including lysosomes. This MAb shows a restricted reactivity to cells of the monocyte/macrophage system. It specifically reacts with blood monocytes and stains resident macrophages in a wide variety of human tissues. This MAb does not stain antigen-presenting cells, e.g., Langerhans cells. Reportedly, its reactivity is restricted to histiocytes and macrophages.

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 0348, Anti-Histiocytoma Marker (D11)

Product attributes

Antibody number	#0348
Antibody reactivity (target)	Histiocytoma Marker
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	D11
Isotype	IgG1, kappa
Molecular weight	125 kDa - 135 kDa
Synonyms	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Membrane preparation from human hepatocytes
Antibody target cellular localization	Cytoplasmic, Plasma membrane
Species reactivity	Human
Antibody application notes	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, Does not react with mouse, rat or pig, others not tested
Positive control	Liver or histiocytoma
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	Immunology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Monocytes/macrophages

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

Rudinskaya TD, Poltoranina VS, Baranov VN, Petrovichev NN, Voytenkov BO, Tretyakov LO. D11, a novel monoclonal antibody specific for human mature macrophages and peripheral blood monocytes. *Immunol Lett.* 1992 Jun;33(1):1-7. | Frenkel MA, Tupitsyn NN, Rudinskaja TD, Poltoranina VS, Lebedeva NB, Volkova MA, Kaletin GI. [Use of a new antimacrophage monoclonal antibody D11 in diagnosis of hemoblastosis]. [Article in Russian] *Gematol Transfuziol.* 1995 Jul-Aug;40(4):13-6. | Tupitsyn NN, et al. Reactivity of anti-macrophage monoclonal antibody D11 in human leukemia and malignant lymphoma. *Int J Cancer.* 1996;68(2):160-3. | Petrovichev NN, et al. Antimacrophage monoclonal antibody D11 in the diagnosis of tumors of histiocytic origin. *Acta Cytol.* 1997;41(2):357-63.?

This datasheet was generated on December 24, 2024 at 05:05:56 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-histiocytoma-marker-d11/>