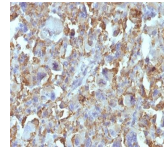


# Histiocytoma Marker Monoclonal Mouse Antibody (D11)



In Western blotting, this antibody detects an antigen of 125 kDa in human liver and 135 kDa in tumors of histiocytic origin.

## 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](mailto: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)**

Antibody #	prefix	Conjugation	Ex/Em	Concentration	Storage Buffer
BNC04	CF®405S		404/431 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC05	CF®405M		408/452 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC88	CF®488A		490/515 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC43	CF®543		541/560 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC55	CF®555		555/565 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC68	CF®568		562/583 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC94	CF®594		593/614 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC40	CF®640R		642/662 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC47	CF®647		650/665 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC61	CF®660R		663/682 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC80	CF®680		681/698 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC81	CF®680R		680/701 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC70	CF®770		770/797 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCR	R-PE (PE)		496, 546, 565/578 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCA	APC		650/660 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCP	PerCP		482/677 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCB	Biotin		N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCAP	Alkaline Phosphatase		N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCH	Horseradish Peroxidase		N/A	0.1 mg/mL	PBS, 0.05% BSA, no azide
BNUB	Purified, with BSA		N/A	0.2 mg/mL	PBS, 0.05% BSA, 0.05% azide
BNUM	Purified, BSA-free		N/A	1 mg/mL	PBS, no BSA, no azide

## 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, Rudinskaya 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.??

Product attributes	
Antibody number	0348
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
Cellular localization	Cytoplasmic, Membrane/cell surface
Species reactivity	Human
Applications	Immunohistology (formalin)
Application notes	Does not react with mouse, rat or pig, others not tested, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
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)
Research areas	Immunology
Cell/tissue expression	Monocytes/macrophages