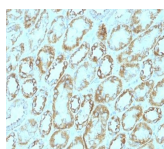


CD171 / L1CAM Monoclonal Mouse Antibody (UJ127)



Product Description

Recognizes a cell surface protein of 220-240 kDa, identified as L1 cell adhesion molecule. The L1CAM gene, which is located in Xq28, is involved in three distinct conditions: 1) HSAS (hydrocephalus-stenosis of the aqueduct of Sylvius); 2) MASA (mental retardation, aphasia, shuffling gait, and adducted thumbs); and 3) SPG1 (spastic paraplegia). The L1, neural cell adhesion molecule (L1CAM) also plays an important role in axon growth, fasciculation, and neural migration as well as in mediating neuronal differentiation. Expression of L1 protein is restricted to tissues arising from neuroectoderm. This MAb is useful in the identification of primitive neuroectodermal tumors. It binds to tumors of neuroectodermal and glial origin e.g. neuroblastoma and Schwannomas. It does not bind to pediatric or adult brain.

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 0679, Anti-CD171 (UJ127)

Product attributes

Antibody number	#0679
Antibody reactivity (target)	CD171, L1CAM
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	UJ127
Isotype	IgG1, kappa
Molecular weight	220-240 kDa
Synonyms	CAML1, HSAS1, Hyd, L1 Cell Adhesion Molecule, L1-NCAM, MASA, MIC5, NCAM-L1, Nerve-growth factor-inducible large external glycoprotein, Neural cell adhesion molecule L1, NILE, S10, SPG1
Human gene symbol	L1CAM
Entrez gene ID	3897
SwissProt	P32004
Unigene	52818
Immunogen	Homogenous suspension of 16 week human fetal brain
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Flow, surface (published for clone), IHC (FFPE) (published for clone), IP (published for clone), WB (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	Neuroblastomas or Schwannomas
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, Cell adhesion, Developmental biology
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	Neuroectodermal 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. Eur J Gynaecol Oncol (2008) 29(1):26-30. (IHC, FFPE; WB)
2. Anticancer Res (2009) 29: 1195-1200. (IHC, FFPE)
3. Mol Biol Cell (2010) 21: 1671-1685. (IHC, FFPE; Flow, IP)

This datasheet was generated on January 28, 2026 at 01:03:34 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd171-l1cam-l1-cell-adhesion-molecule-uj127-11/>