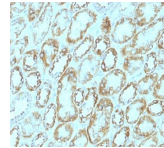


CD171 / L1CAM Monoclonal Mouse Antibody (UJ127)



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)**

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)

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 |
| 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 |
| 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, LI-COR Bioscience.