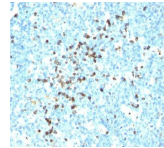


# Human Lambda Light Chain Monoclonal Mouse Antibody (LcN-2)



## Product Description

This MAb is specific to lambda light chain of immunoglobulin and shows no cross-reaction with lambda light chain or any of the five heavy chains. In mammals, the two light chains in an antibody are always identical, with only one type of light chain, kappa or lambda. The ratio of Kappa to Lambda is 70:30. However, with the occurrence of multiple myeloma or other B-cell malignancies this ratio is disturbed. Antibody to the lambda light chain is reportedly useful in the identification of leukemias, plasmacytomas, and certain non-Hodgkin's lymphomas. Demonstration of clonality in lymphoid infiltrates indicates that the infiltrate is malignant. 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.

## Product attributes

|  |  |
|--|--|
| <b>Antibody number</b>                       | #0631  |
| <b>Antibody reactivity (target)</b>          | Human Lambda Light Chain   |
| <b>Antibody type</b>                         | Anti-Human Immunoglobulin, Primary   |
| <b>Host species</b>                          | Mouse  |
| <b>Clonality</b>                             | Monoclonal   |
| <b>Clone</b>                                 | LcN-2  |
| <b>Isotype</b>                               | IgG2a, kappa   |
| <b>Molecular weight</b>                      | ~22.5 kDa  |
| <b>Synonyms</b>                              | Bence Jones Protein; BJP; IGLC 1/2/3; Mcg Marker; Paraprotein  |
| <b>Human gene symbol</b>                     | IGL  |
| <b>Entrez gene ID</b>                        | 3535 & 3546  |
| <b>SwissProt</b>                             | P01701 & P01842  |
| <b>Unigene</b>                               | 449585   |
| <b>Immunogen</b>                             | Purified human IgG   |
| <b>Antibody target cellular localization</b> | Plasma membrane, Secreted (extracellular)  |
| <b>Verified antibody applications</b>        | IF (verified), IHC (FFPE) (verified), WB (verified)  |
| <b>Species reactivity</b>                    | Human  |
| <b>Antibody application notes</b>            | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.25-0.5 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, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user |
| <b>Positive control</b>                      | 293T, Raji or hPBL cells. Tonsil or Spleen   |
| <b>Shipping condition</b>                    | Room temperature   |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended  |
| <b>Regulatory status</b>                     | For research use only (RUO)  |
| <b>Antibody/conjugate formulation</b>        | 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  |
| <b>Cell/tissue expression</b>                | B-cells  |
| <b>Antibody research areas</b>               | Cancer, Immunology, Inflammation   |
| <b>Tumor expression</b>                      | Leukemia/lymphoma  |



## Catalog Number Key for Primary Antibody Conjugates

| Antibody # prefix | Conjugation           | Ex/Em (nm) | Laser line | Detection channel        | Dye Features                     |
|-------------------|-----------------------|------------|------------|--------------------------|----------------------------------|
| BNC04             | CF@405S               | 404/431    | 405        | DAPI (microscopy), AF405 | <a href="#">CF@405S Features</a> |
| BNC88             | CF@488A               | 490/515    | 488        | GFP, FITC                | <a href="#">CF@488A Features</a> |
| BNC68             | CF@568                | 562/583    | 532, 561   | RFP, TRITC               | <a href="#">CF@568 Features</a>  |
| BNC94             | CF@594                | 593/614    | 561        | Texas Red®               | <a href="#">CF@594 Features</a>  |
| BNC40             | CF@640R               | 642/662    | 633-640    | Cy@5                     | <a href="#">CF@640R Features</a> |
| BNC47             | CF@647                | 650/665    | 633-640    | Cy@5                     | <a href="#">CF@647 Features</a>  |
| BNCB              | Biotin                | N/A        | N/A        | N/A                      |                                  |
| BNUB              | Purified              | N/A        | N/A        | N/A                      |                                  |
| BNUM              | Purified,<br>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 GE Healthcare; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

---

This datasheet was generated on September 30, 2022 at 06:02:22 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/monoclonal-anti-lambda-light-chain-lcn-2/>