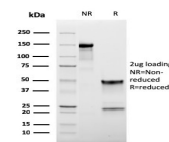


# CD62L / L-Selectin Monoclonal Mouse Antibody (LAM1-116)



## Product Description

Selectins, also designated CD62 antigens, comprise a family of carbohydrate-binding proteins involved in mediating cellular interactions with leukocytes. L-Selectin (also designated LECAM-1 or CD62L) is expressed on the majority of B and naive T cells and on most monocytes, neutrophils and eosinophils. L-Selectin interacts with specific carbohydrates expressed by activated endothelial cells. P-Selectin (also designated GMP-140 or CD62P), expressed on activated platelets and endothelial cells, and E-Selectin (also designated ELMA-1 or CD62E), expressed on endothelial cells, exhibit overlapping ligand specificities. Both recognize sialyl-Le (x) as a ligand and bind to specific carbohydrates on neutrophils and monocytes.

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 1587, Anti-CD62L|L-Selectin (LAM1-116)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

|  |  |
|--|--|
| <b>Antibody number</b>                       | #1587  |
| <b>Antibody reactivity (target)</b>          | CD62L, L-Selectin  |
| <b>Antibody type</b>                         | Primary  |
| <b>Host species</b>                          | Mouse  |
| <b>Clonality</b>                             | Monoclonal   |
| <b>Clone</b>                                 | LAM1-116   |
| <b>Isotype</b>                               | IgG1, kappa  |
| <b>Molecular weight of antigen</b>           | 65-75 kDa  |
| <b>Synonyms</b>                              | A.11; AI528707; CD62 antigen ligand; CD62L; gp90-MEL; IgA nephropathy; susceptibility to; included; L Selectin; L-selectin; LAM1; LECAM1; LEU8; Leukocyte surface antigen Leu-8; Leukocyte-endothelial cell adhesion molecule 1; Lnhr; LSEL; Ly-22; LYAM1; Lymph node homing receptor; Lymphocyte antigen 22; Lymphocyte surface MEL-14 antigen; Pln homing receptor; PLNHR; SELL; TQ1 |
| <b>Human gene symbol</b>                     | SELL   |
| <b>Entrez gene ID</b>                        | 6402   |
| <b>SwissProt</b>                             | P14151   |
| <b>Unigene</b>                               | 728756   |
| <b>Immunogen</b>                             | MAb raised against supernatant from phorbol myristic acid activated human peripheral blood leukocytes.   |
| <b>Antibody target cellular localization</b> | Plasma membrane  |
| <b>Species reactivity</b>                    | Human  |
| <b>Expected antibody applications</b>        | ELISA (published for clone), Flow, surface (published for clone), Functional studies (published for clone), IHC (FFPE) (published for clone), IP (published for clone), WB (published for clone)   |
| <b>Antibody application notes</b>            | ELISA: For coating, order antibody without BSA; Flow cytometry: 0.5-1 ug/million cells; Western blot 0.5-1.0 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody   |
| <b>Positive control</b>                      | HPBL or Jurkat whole cell lysate.  |
| <b>Shipping condition</b>                    | Room temperature   |
| <b>Storage Conditions</b>                    | Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light  |
| <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>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended  |
| <b>Product origin</b>                        | 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.  |
| <b>Cell/tissue expression</b>                | Eosinophils, Lymphocytes, Monocytes/macrophages, Neutrophils   |

| 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>  |
| BNC74             | CF@740                | 742/767    | 633-685    | 775/50                   | <a href="#">CF@740 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 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. J Clin Invest (1997) 99(8): 2036-2044. (Flow)
2. J Clin Invest (1997) 100(5): 996-1005. (functional studies)
3. J Biol Chem (1997) 272(21): 13961-13965. (Flow, IP)
4. Ann Rheum Dis (2000) 59:297-299. (ELISA)
5. J Immunol (2002) 169:2034-2043. (ELISA)
6. Immunity (2003) 19(5): 713-724. (WB)
7. Fert Stert (2007) 88(4) Supplement: 1207-1211. (IHC, FFPE)
8. J Leukocyte Biol (2008) 83: 99-105. (Flow, IF)

This datasheet was generated on June 17, 2026 at 03:00:55 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd62l-l-selectin-monoclonal-mouse-antibody-lam1-116/>