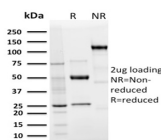


Langerin / CD207 Recombinant Monoclonal Rabbit Antibody (LGRN/3136R)



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

Product Description

Langerhans cells (LCs) are a subset of immature dendritic cells (DCs) that specifically localize in the epidermis and other mucosal epithelia. Epidermal LCs possess strong immuno-stimulatory capacity and play a central role in the initiation and regulation of immune responses. Langerin (CD207) is a Ca²⁺-dependent, C-type lectin domain containing, type II transmembrane protein that induces epidermal LCs to differentiate into Birbeck granules (BG). BGs are organelles with superimposing and zipper membranes that influence proper class I type antigen presentation to the circulating T cells. Human spleen, lymph node, thymus, liver, lung and heart express Langerin protein.

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 3136, Anti-CD207|Langerin (LGRN/3136R)

Product attributes

| | |
|--|---|
| Antibody number | #3136 |
| Antibody reactivity (target) | CD207, Langerin |
| Antibody type | Primary |
| Host species | Rabbit |
| Clonality | Monoclonal, Recombinant Monoclonal |
| Clone | LGRN/3136R |
| Isotype | IgG, kappa |
| Molecular weight of antigen | 48 kDa |
| Synonyms | C-type lectin domain family 4 member K; CD207; CLEC4K; Langerhans cell specific c type lectin; Langerin |
| Human gene symbol | CD207 |
| Entrez gene ID | 50489 |
| SwissProt | Q9UJ71 |
| Unigene | 199731 |
| Immunogen | Recombinant fragment (around aa74-213) of human Langerin protein (exact sequence is proprietary) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Antibody application notes | For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |
| Positive control | HeLa or A549 cells. Skin. |
| 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 |
| 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 |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Product origin | Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Cell/tissue expression | Heart, Lymph node, Dendritic cells, Liver, Lung, Spleen, Thymus |

| 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.

This datasheet was generated on June 15, 2026 at 01:31:02 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/langerin-cd207-recombinant-monoclonal-rabbit-antibody-lgrn-3136r/>