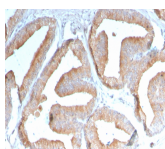


ICOS-L / ICOS Ligand / B7RP-1 Monoclonal Mouse Antibody (ICOSL/3111)



Product Description

This MAbs recognizes a protein of 36 kDa, identified as ICOS-L. It is a ligand for the ICOS receptor that initiates T and B cell proliferation and cytokine secretion. ICOS-L interactions play an essential role in T cell-dependent B cell activation in peripheral lymphoid organs such as spleen and lymph nodes. ICOS-L protein is present in myeloid leukocytes, and by Northern blot there are 2.4, 3.0, and 7.0 kb transcripts in brain, heart, kidney, and liver, with lower expression in colon and thymus, and a 1.1 kb transcript in leukocytes. Tumor necrosis factor alpha (TNF), granulocyte-macrophage colony-stimulating factor (GM-CSF) and interleukin-4 (IL-4) enhance its expression. LPS-induced up-regulation of ICOS-L is dependent on the MyD88-dependent signaling pathway.

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 3111, Anti-B7RP-1|ICOS Ligand|ICOS-L (ICOSL/3111)

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

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #3111 |
| Antibody reactivity (target) | B7RP-1, ICOS Ligand, ICOS-L |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | ICOSL/3111 |
| Isotype | IgG1, kappa |
| Molecular weight | 36 kDa |
| Synonyms | B7 homolog 2 (B7H2); B7 like protein GI50; B7 related protein 1 (B7RP1); CD275; CD275 antigen; GL50; ICOS LG; ICOS ligand (ICOSL); Inducible T cell co stimulator ligand; LICOS; Transmembrane protein B7 H2 ICOS ligand |
| Human gene symbol | ICOSLG |
| Entrez gene ID | 23308 |
| SwissProt | O75144 |
| Unigene | 14155 |
| Immunogen | Recombinant fragment of human ICOSLG (B7RP-1) protein (around aa 23-149) (exact sequence is proprietary) |
| Verified antibody applications | Flow (verified), IF (verified), IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | U937, HL60, HuT78 and HUVEC lysates (WB). Human tonsil tissue (IHC). |
| 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 |
| Validated in protein array | Monospecific |
| 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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Cell/tissue expression | Myeloid cells |

| 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 January 14, 2026 at 09:57:17 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/icos-l-icos-ligand-b7rp-1-monoclonal-mouse-antibody-icosl-3111/>