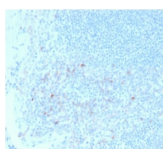


## CD25/ IL2RA Monoclonal Mouse Antibody (IL2RA/2395)



### Product Description

This antibody recognizes a protein of 55 kDa, identified as CD25. It is expressed on activated T- and B-cells and activated monocytes/macrophages. With respect to lymphomas, CD25 is present on malignant cells of Hodgkin's disease, HTLV-1 adult T-cell leukemia, cutaneous T-cell lymphoma, and hair cell leukemia. Increased levels of soluble CD25 are observed in the leukemias/lymphomas and inflammatory/ autoimmune diseases. CD25 molecule alone appears to function as a low affinity receptor and associates with CD122 (IL-2R  $\alpha$  chain, p75) and CD132 (common chain) to form the high affinity IL-2 receptor complex. CD25 antibodies detect three epitope regions, A, B and C. This MAAb This antibody recognizes the epitope B, which is located at residue 3-104 of CD25 and does not block IL-2 binding to CD25.

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 2395, Anti-CD25|IL2RA (IL2RA/2395)**

### Product attributes

Antibody number	#2395
Antibody reactivity (target)	CD25, IL2RA
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IL2RA/2395
Isotype	IgG2b
Molecular weight of antigen	55 kDa
<b>Synonyms</b>	IL2 Receptor alpha; IL2R alpha chain; IL2R; IL2R subunit alpha IL2RA; Interleukin-2 receptor subunit alpha; Ly43; p55; p55 chain; T Cell Growth Factor Receptor; TAC antigen; TCGFR
Human gene symbol	IL2RA
Entrez gene ID	3559
SwissProt	P01589
UniGene	231367
Immunogen	Recombinant fragment of human IL2RA protein (around aa 42-183) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 2-4 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user
Positive control	PHA-stimulated human lymphocytes. Human lymph node or tonsil
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
Validated in protein array	Monospecific
Antibody research areas	Immunology
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	B-cells, Monocytes/macrophages, T-cells
Tumor expression	Leukemia/lymphoma

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, 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 6, 2026 at 09:36:37 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd25-il2ra-monoclonal-mouse-antibody-il2ra-2395/>