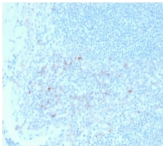


# CD25/ IL2RA Monoclonal Mouse Antibody (IL2RA/2395)



## 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  $\alpha$ - $\beta$  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 MAb 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.

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	55 kDa
Synonyms	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 assay	Monospecific
Antibody research areas	Immunology
Cell/tissue expression	B-cells, Monocytes/macrophages, T-cells
Tumor expression	Leukemia/lymphoma