

IL-2R α (CD25) Recombinant monoclonal antibody (Rabbit IgG κ)

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Interleukin-2 receptor subunit alpha, IL-2 receptor subunit alpha, IL-2R subunit alpha, IL-2-RA, TAC antigen, p55
Application	Flow Cytometry, FUNC, IF, IHC, WB
Clone	Basiliximab
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	Basiliximab was prepared by immunizing BALB/c mice with CTC 16 cells.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Recognizes human IL-2R and binds to the epitope E ¹¹⁶ RIYHFV ¹²² within the extracellular domain of the CD25 subunit. Binding is abolished when 2 or more residues within the epitope are mutated. This epitope overlaps with the interaction site of CD25 and IL-2. CD25 is found in the high-affinity IL-2R ($\alpha\beta\gamma$) – T cells only express CD25 upon activation. IL-2 binding to IL-2R on T cells modulates proliferation and differentiation.

Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original Human IgG1 format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P01589

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com