IL-2Rα (CD25) Recombinant monoclonal antibody (Rabbit IgGκ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS546-0200

200µg

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

ImmunogenBasiliximab 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 E¹¹⁶

RIYHFV¹²² 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 ($\alpha\beta\gamma c$) – 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



info-

eu@enzolifesciences.com