

# IL-2R $\alpha$ (CD25) Recombinant monoclonal antibody (Rabbit IgG $\kappa$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS546-0200	200 $\mu$ g
-----------------	-------------

## 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 <sup>116</sup> RIYHFV <sup>122</sup> 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 ( $\alpha\beta\gamma\epsilon$ ) – T cells only express CD25 upon activation. IL-2 binding to IL-2R on T cells modulates proliferation and differentiation.

<b>Technical Info / Product Notes</b>	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.
<b>UniProt ID</b>	P01589
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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