

# IL-2R $\alpha$ (CD25) Recombinant monoclonal antibody (Human IgG1 $\kappa$ )

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Interleukin-2 receptor subunit alpha, IL-2 receptor subunit alpha, IL-2R subunit alpha, IL-2-RA, TAC antigen, p55
<b>Application</b>	Flow Cytometry, FUNC
<b>Clone</b>	Daclizumab
<b>Formulation</b>	Liquid. In PBS containing 0.02% Proclin 300.
<b>Host</b>	Human
<b>Immunogen</b>	Cultured T-cells derived from a patient with Mycosis fungoides.
<b>Isotype</b>	IgG1κ
<b>Purity Detail</b>	Protein A affinity purified.
<b>Species Reactivity</b>	Human
<b>Specificity</b>	Recognizes the p55 chain of IL2R on human T cells, Tac antigen, which is expressed on T cells activated by mitogens, soluble antigens, and alloantigens but not on resting T cells, thymocytes, B cells, monocytes, activated B cells, leukemic T cell blasts from patients with acute lymphoblastic leukemia, and long-term T cell lines.
<b>UniProt ID</b>	P01589

Last modified: May 29, 2024



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)