

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

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

## Ordering Information

[Order Online »](#)

ENZ-ABS545-0200	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
Clone	Daclizumab
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	Cultured T-cells derived from a patient with Mycosis fungoides.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	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.
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](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)