

V α TCR Recombinant monoclonal antibody (34) (Mouse IgG1 κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS606-0200	200 μ 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	TCR variable α chain, T-cell receptor variable α chain
Application	ELISA, Flow Cytometry, WB
Clone	34
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Purified recombinant soluble 85.33 scTCR.
Isotype	IgG1k
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Technical Info / Product Notes	This full-length, reformatted mouse antibody was made using the variable domain sequences of the original Mouse scFv format, for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P01848



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

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

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

France
Phone: +33 1 402 44 555
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

This antibody is covered by our **Worry Free Guarantee**