

# Integrin $\alpha 4$ Recombinant monoclonal antibody (Rabbit IgG $\kappa$ )

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

## Ordering Information

[Order Online »](#)

ENZ-ABS642-0200	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	CD49d, α4 subunit of VLA-4 receptor, ITGA4, Integrin α-IV, Natalizumab
Application	Flow Cytometry, FUNC
Clone	Natalizumab
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	The RAMOS cell line, a non-EBV releasing cell line, injected into mice.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Recognizes the α4 subunit of α4β1 integrin (also known as very late antigen 4 [VLA-4] or CD49d-CD29) and α4β7 integrin.
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original human IgG4 format, for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P13612

Worry-free Guarantee

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



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)