## Integrin α4 Recombinant monoclonal antibody (Rabbit IgGκ)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS642-0200

200µg

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** IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

**Specificity** Recognizes the  $\alpha 4$  subunit of  $\alpha 4\beta 1$  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.

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com