

# CD51 (human) monoclonal antibody (L230)

CD51 (integrin  $\alpha$ V subunit) associates with CD61 (integrin  $\beta$ 3 subunit) binding to the RGD sequence in extracellular matrix proteins like vitronectin, fibrinogen, thrombospondin, von Willebrand factor and osteopontin. The  $\alpha$ V/ $\beta$ 3 heterodimer is expressed in endothelial cells, leukocytes, NK cells, macrophages, neutrophils, platelets and several malignant tumor cells and seems to play a role in embryo implantation, angiogenesis, wound healing and cancer.  $\alpha$ V/ $\beta$ 3 is also highly expressed in osteoclasts being important for bone resorption.

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

Citations: 15

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |             |
|------------------|-------------|
| ALX-803-304-C100 | 100 $\mu$ g |
|------------------|-------------|

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

|                    |   |
|--------------------|---|
| Handling           | Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C. |
| Short Term Storage | +4°C  |
| Long Term Storage  | -20°C   |
| Shipping           | Blue Ice  |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                      |  |
|----------------------|--|
| Alternative Name     | αV integrin, Vitronectin receptor α subunit                            |
| Application          | ELISA, Flow Cytometry, FUNC, IF, IHC, IP                               |
| Application Notes    | Functional Application: αV integrin blocking                           |
| Clone                | L230   |
| Crossreactivity      | Does not cross-react with other species.                               |
| Formulation          | Liquid. In PBS containing 0.05% sodium azide.                          |
| Host                 | Mouse  |
| Immunogen            | Human endothelial cells.   |
| Isotype              | IgG1   |
| Purity Detail        | Protein A-affinity purified.   |
| Species Reactivity   | Human  |
| Specificity          | Recognizes heterodimers αVβ3, αVβ1, αVβ5 and αVβ6.                     |
| UniProt ID           | P06756   |
| Worry-free Guarantee | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> . |



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)