

# CD51 polyclonal antibody

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: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-537-R100	100 $\mu$ l
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Vitronectin receptor  $\alpha$  subunit,  $\alpha$ V integrin

**Application** ELISA, ICC, IHC (FS), IHC (PS), IP, WB

**Formulation** Liquid. Neat serum containing 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to C-terminal portion of human CD51 ( $\alpha$ V integrin) (aa 1021-1034) (cytoplasmic domain).

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:1,000)Immunohistochemistry (frozen sections, 1:500)Immunohistochemistry (paraffin sections, 1:500)Immunoprecipitation (5 $\mu$ l per 5 $\times$ 10<sup>6</sup> cells)Western Blot (1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Species independent

**Specificity** Recognizes the cytoplasmic domain of CD51.

**UniProt ID** P06756

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