

CD51 polyclonal antibody

CD51 (integrin αV 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 »](#)

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 α subunit, α 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 (α 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 μ l per 5 \times 10⁶ 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](#)