

CD51 polyclonal antibody

CD51 (integrin α V subunit) associates with CD61 (integrin β 3 subunit) binding to the RGD sequence in extracellular matrix proteins like vitronectin, fibrinogen, thrombospondin, von Willebrand factor and osteopontin. The α V/ β 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. α V/ β 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 μ 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 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 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.

Dilutions/Conditions

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](#).



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)