

Vitronectin monoclonal antibody (A27)

Vitronectin is a plasma glycoprotein that circulates in the blood as a mixture of both 75kDa and 65kDa forms. Vitronectin is a major cell adhesive glycoprotein, and a common component of extracellular matrix and plasma. It competes effectively with other plasma proteins, and often participates in cell attachment, regulation of blood coagulation and immune responses. It has similar tissue distribution to fibronectin, which is recognized by its integrin receptor.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Serum spreading factor, S-protein
Application	Affinity chromatography, ELISA, IHC (FS), IHC (PS), WB
Clone	A27
Crossreactivity	Does not cross-react with human or horse. There is no evidence for cross-reactivity with other connective tissue proteins (fibronectin, elastin, collagen, laminin).
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	507525
Host	Mouse
Immunogen	Lysed bovine corneal endothelial cells and extracellular matrix.
Isotype	IgG1k
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Bovine, Sheep

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024