

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](#).

Ordering Information

[Order Online »](#)

BPD-CSI-004-27-1	1mg
BPD-CSI-004-27-04	400µg

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 IgG1κ

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-be@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-
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)