

Fibronectin (bovine) monoclonal antibody (A22)

Fibronectin is an adhesive glycoprotein with a molecular mass of 440kDa which appears important in the formation of a provisional matrix to promote cell adhesion and migration during wound healing. Its age-dependent increase in plasma and tissues may accompany a proteolytic breakdown by a number of neutral proteases in pathological states, especially in tumor growth. Several of its proteolytic breakdown products exhibit unexpected and primarily harmful biological activities.

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

Ordering Information

[Order Online »](#)

BPD-CSI-005-22-1	1mg
BPD-CSI-005-22-04	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 FN

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

Clone A22

Crossreactivity Does not cross-react with human. There is no evidence for cross-reactivity with other connective tissue proteins (vitronectin, elastin, collagen, laminin).

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

GenBank ID 280794

Host Mouse

Immunogen Bovine corneal endothelial cells on beads.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Bovine

UniProt ID P07589

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-
be@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)