

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-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)