

# Fibronectin monoclonal antibody (A17)

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-005-17-1	1mg
BPD-CSI-005-17-04	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
Application Notes	The antibody inhibits integrin-mediated cell adhesion to the cell binding domain of fibronectin.
Clone	A17
Crossreactivity	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	Lysed bovine corneal endothelial cells and extracellular matrix.
Isotype	IgG1κ
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Bovine, Chicken, Human
Specificity	Epitope is located in the 120kD cell binding fragment.
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)