

# Fibronectin monoclonal antibody (A32)

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](#).

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	FN
<b>Application</b>	ELISA, IHC (FS), IHC (PS), IP, WB
<b>Application Notes</b>	The antibody inhibits cell adhesion to fibronectin mediated by the heparin-binding domain.
<b>Clone</b>	A32
<b>Crossreactivity</b>	There is no evidence for cross-reactivity with other connective tissue proteins (vitronectin, elastin, collagen, laminin).
<b>Formulation</b>	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
<b>GenBank ID</b>	280794
<b>Host</b>	Mouse
<b>Immunogen</b>	Bovine corneal endothelial cells.
<b>Isotype</b>	IgG1k
<b>Purity Detail</b>	Protein A/G-purified.
<b>Source</b>	From culture supernatant.
<b>Species Reactivity</b>	Bovine, Human
<b>UniProt ID</b>	P07589

Last modified: May 29, 2024



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)