

Fibronectin (plasma) (human) monoclonal antibody (BF12)

Fibronectins are high molecular weight, disulphide-linked, dimeric cell adhesion glycoproteins found in basement membranes and in the interstitial connective tissue matrix. A single fibronectin gene is subject to alternative splicing in a cell-type-, development- and age-regulated manner which gives rise to multiple molecular forms. In addition to their prominent role in adhesion, fibronectins have been reported to mediate various aspects of cellular interaction, including migration during development and wound-healing, haemostasis, and the regulation of cell growth and differentiation. Cellular fibronectins (cFn) are found in low amounts in normal human plasma and tissues, but they are abundant in the plasma of carcinoma patients and in the stroma of various carcinomas. In contrast, a soluble form of fibronectin produced by hepatocytes is readily detectable in plasma and becomes deposited in pericellular matrices and within tissues. This form of fibronectin, referred to as 'plasma fibronectin' (pFn), differs from cFn by the absence of an amino acid sequence, known as extra domain A1.

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

Ordering Information

[Order Online »](#)

BML-FG6020-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name pFn

Application ICC, WB

Application Notes Detects a band of ~225-250kDa by Western blot.

Clone BF12

Formulation Liquid. In PBS containing 0.1% sodium azide and 1mg/mL BSA.

Host Mouse

Immunogen Purified fibronectin from A8387 fibrosarcoma cells.

Isotype IgG1

Purity Detail Protein A affinity chromatography.

Species Reactivity Human

Specificity Recognizes the cell binding region of fibronectin.

UniProt ID P02751

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

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-
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)