

Fibronectin

From human plasma
Lyophilizate, before lyophilization filtered through a 0.2 µm pore size membrane.

Cat. No. 10 838 039 001 5 mg

Version 22
Content version: October 2018

Store at +2 to +8°C

Product overview

Formulation Lyophilized (contains glycine and sodium chloride); before lyophilization filtered through a 0.2 µm pore size membrane.

Preparation Human fibronectin is purified from human plasma (1). The raw material used for the production is negative as tested for Anti-HIV antibodies and HBs-antigen.

Primary structure Human plasma fibronectin consists of two similar polypeptide chains (α-, β-chain, each of 220 kDa), connected by two disulfide bridges (2-4).

Properties

Molecular weight	440,000 Da
Purity	> 95% pure as determined by HPLC, additionally checked on SDS-PAGE.
Species specificity	effective on most mammalian cells.
Biological activity	Tested for the promotion of adherence and growth of human umbilical vein endothelial cells.

Description

Fibronectin is a large dimeric glycoprotein. Fibronectin is widely distributed in a soluble form (plasma fibronectin) in plasma and other body fluids. Many cell types synthesize and secrete fibronectin, but most circulating fibronectin is produced by hepatocytes. Fibronectin is also widely distributed in an insoluble form in tissues (cellular fibronectin), where it is covalently cross-linked into multimeric fibers (2). Plasma fibronectin is not identical to that in extracellular matrices and on cell surfaces (cellular fibronectin), but is equally active in cell attachment. Fibronectin has several adhesive functions, e.g., cell-to-cell attachment, cell adherence to plastic or basement membranes and clot stabilization. Each polypeptide chain of fibronectin can be divided into functional domains (e.g., cell binding domain, collagen binding domain, heparin binding domain) (2-4).

Application

Fibronectin promotes the attachment and subsequent spreading of many cells. The cell adhesion region (cell binding domain) interacts with mammalian cells and promotes their binding to and spreading on plastic. Fibronectin also binds to other extracellular and basement membrane components and mediates cell attachment to collagen (5-10).

Reconstitution

Add 5 ml sterile water to yield a final concentration of 1 mg/ml. Incubate for 30 - 60 min at +37°C to dissolve. **Do not agitate!** Upon reconstitution, the solution may contain a small amount of insoluble aggregated material. Occurrence of these aggregates is a phenomenon inherent to fibronectin (1) and does not influence product performance.

Working concentration

For the coating of cell culture vessels 5 µg/cm² is used.

Storage/Stability

The lyophilizate is stable at +2 to +8°C, the reconstituted solution is stable at -15 to -25°C.

Note: It is recommended to store the reconstituted solution in aliquots; at -15 to -25°C. Repeated freezing and thawing should be avoided.

Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

Coating cell culture vessels with fibronectin

Protocol

Please refer to the following table.

Step	Action
1	Dissolve the lyophilizate with sterile tissue grade water to a final concentration of 1 mg/ml as described above (reconstitution).
2	Dilute an appropriate aliquot with sterile PBS or basal medium to a final concentration of 50 µg/ml.
3	Pipette 100 µl of this solution (50 µg/ml) per 1 cm ² surface area to be coated (5 µg/cm ²). Note: This can be increased or decreased to fit the application. However, the solution should completely wet the surface.
4	Incubate for about 45 min at +15 to +25°C.
5	Remove the solution carefully from the edges without touching the surface area with the pipette.
6	It is possible, but not necessary, to wash the coated surface with medium or buffer. Note: The plates should not be allowed to dry! The vessels should be used immediately.

References

- 1 Vuento, M. & Vaheri, A. (1979) *Biochem. J.* **183**, 331-337.
- 2 Proctor, R. A. (1987) *Rev. Inf. Dis.* **9**, 317-321.
- 3 Mosher, D. F. (1984) *Ann. Rev. Med.* **35**, 561-575.
- 4 Hynes, R. (1985) *Ann. Rev. Cell Biol.* **1**, 67-90.
- 5 Grinnel, F. (1981) In: *Tissue Growth Factors* (Baserga, R., ed.) Springer Verlag, Berlin, Heidelberg, New York, pp. 277-299.
- 6 Ruoslahti, E. et al. (1982) *Methods Enzymol.* **82**, 803-831.
- 7 Kleinman, H. K., Klebe, R. J. & Martin, G. R. (1981) *J. Cell Biol.* **88**, 473-485.
- 8 Yamada, K. M. (1983) *Ann. Rev. Biochem.* **52**, 761-799.
- 9 Kleinman, H. K. et al. (1987) *Anal. Biochem.* **166**, 1-13.
- 10 McKeown-Longo, P. J. (1987) *Rev. Inf. Dis.* **9**, 322-334.

Changes to Previous Version

Editorial changes.

Trademarks

All third party product names and trademarks are the property of their respective owners.

Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit sigma-aldrich.com, and select your home country. Country-specific contact information will be displayed.