

Cell Attachment Factors

PromoCell

Instruction Manual

Product	Size	Catalog Number
Fibronectin Solution, bovine (1 mg/ml)	5 ml	C-43050
Fibronectin Solution, human (1 mg/ml)	5 ml	C-43060

Description

Fibronectin is an attachment factor that facilitates the attachment and cytoplasmic spreading of all types of anchorage-dependent cells. Fibronectin is particularly useful for cells that are not capable of synthesizing their own biomatrix or when culturing cells in serum-free medium.

PromoCell Fibronectin is a clear sterile solution obtained by affinity purification on gelatin-sepharose from irradiated citrated bovine or human plasma, respectively. The Fibronectin solutions contain 1 mg/ml Fibronectin and buffer salts.

Instructions for Use

Depending on the cell type and culture conditions, dilute the Fibronectin So-

lution to 1 - 10 µg/ml final concentration in Dulbecco's PBS w/o Calcium and Magnesium (C-40232). Overlay the culture surface of your tissue culture vessel with an amount of the diluted Fibronectin Solution sufficient to effectively coat the complete surface. Be sure that the entire surface is covered. Place flasks on a level surface at RT for 60 min. Aspirate the excess Fibronectin Solution and use immediately or let air-dry the open vessel under a laminar flow bench. Unused vessels may be stored at 4°C for up to 2 weeks.

Storage and Stability

Store at 4 to 8°C in the dark immediately after arrival. Do not freeze. If stored properly, the product is stable until the expiry date stated on the label.

Quality Control

All lots of PromoCell cell attachment factors are subjected to comprehensive quality control tests. Each lot is routinely tested for coating efficiency, absence of cytotoxicity and promotion of cell attachment and growth. Approved in-house lots are used as a reference.

In addition, all lots have been tested for the absence of microbial contaminants (fungi, bacteria).

Intended Use

The products are for *in vitro* research use only and not for diagnostic or therapeutic procedures. For safety precautions please see appropriate MSDS.