

cFN

native, cellular Fibronectin
human, foreskin fibroblasts

Cat. No.	Amount
PR-389	250 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in Phosphate buffer pH 7.5 and NaCl.

Application

- 1) structural studies: amino acid composition and sequence; folding, configuration
- 2) cell culture: cell attachment, growth, relation to extracellular matrix, regulation of cell migration
- 3) animal studies: dynamics of wound healing, burns, eye abrasions; physiology of metastasis; controls of embryologic development
- 4) mechanism of collagen contraction.

Purity

> 95% by SDS-PAGE

Description

Cellular Fibronectin is different from the usual plasma FN. The cFN molecule contains some domains not present in pFN. Plasma FN is made by liver cells and released into the blood, where it takes part in clotting. Cellular FN is made by general cells of the body, mostly fibroblasts in skin and elsewhere, and becomes part of the extracellular matrix. Plasma FN is involved in early response to a wound (clot formation), while cFN takes part in later stages of wound healing. Malignant cells respond differently to pFN and cFN, and collagen gel contraction is mediated by cFN and not by pFN.

Purified from human foreskin fibroblasts grown in culture.

Selected References:

- Kuschel *et al.* (2006) Cell adhesion profiling using extracellular matrix protein microarrays. *Biotechniques*. **40**:523.
- Zand *et al.* (2003) Differential effects of cellular fibronectin and plasma fibronectin on ovarian cancer cell adhesion, migration, and invasion. *In Vitro Cell Dev Biol Anim.* **39**:178.
- Lee *et al.* (2001) Comparing transforming growth factor beta-2 and fibronectin as pleurodesing agents. *Respirology*. **6**:281.
- Rodrigues *et al.* (2001) Conformational regulation of the fibronectin binding and $\alpha 3 \beta 1$ integrin-mediated adhesive activities of thrombospondin-1. *J Biol Chem.* **276**:27913.
- Yoshizato *et al.* (1999) The interaction of cellular fibronectin with collagen during fibroblast-mediated contraction of collagen gels. *J Invest Dermatol Symp Proc.* **4**:190.
- Jarnagin *et al.* (1994) Expression of variant fibronectins in wound healing: cellular source and biological activity of the EIIIA segment in rat hepatic fibrogenesis. *J Cell Biol.* **127**:2037.