



GIBCOBRL



LIFE TECHNOLOGIES™

Fibronectin (Murine)

Cat. No.: 12173-019

Size: 1 mg

Store at -70°C.

Note: Do not store in dilute solution. Do not freeze/thaw repeatedly.

Form: Liquid

Description:

This preparation was purified from mouse plasma by affinity chromatography on gelatin-Sepharose.

It is supplied at a concentration of 1 mg/ml, determined by light absorption at 280 nm, using $A_{280} = 1.3$ of a 1 mg/ml solution.

Fibronectin has a molecular weight of 220 kDa (reduced) by SDS-PAGE.

This fibronectin is suitable for use in quantitative and qualitative immunochemical methods such as ELISA, Ouchterlony, and immunoblotting.

Doc Rev.: 01/2901

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE™ (800) 828-6686.

Directions for Use:

Mouse fibronectin is supplied frozen in a 2.7 mM potassium chloride, 340 mM sodium chloride, 8 mM sodium phosphate, 1.5 mM potassium phosphate, pH 7.3, and 10% glycerol buffer, with no preservatives added.

Avoid repeated freeze/thaw cycles.

Quality Control:

Determined to be $\geq 95\%$ pure by SDS-PAGE.

References:

Engvall, E. and Ruoslahti, E. (1977) *Int. J. Cancer*, **20**, 1.

Ruoslahti, E. *et al.*, (1981) *Biochem. J.*, **193**, 295.