



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Human Fibronectin 1 Recombinant

Catalog #: 90143

Lot#: 170731

Description: Recombinant human FN is a major component of the provisional ECM produced during sprouting angiogenesis. FN, also known as CD36, is a disulfide-linked monomeric protein consisting of 220 amino acids and migrates as an approximately 28 kDa protein under reducing and non-reducing conditions.

Source: Optimized DNA sequence encoding fibronectin (LYS2175 - GLU2386) including a C-terminal His tag was expressed in *HEK293 cells*.

Formulation: Lyophilized from from 0.2 μ m filtered PBS solution, pH7.2 , 5% Trehalose. at 0.1-1.0 mg/mL in distilled water. This solution can then be diluted into other buffers. To maximize product collection from vial surface, vortex briefly and then spin down to recollect the liquid.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >95%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/ μ g (1EU/ μ g), using the LAL gel clot method.

References:

1. *J. Biol. Chem.*, Feb 2014; 289: 4941 - 4951.
2. *Carcinogenesis*, Jan 2014; 35: 184 - 191.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com