

Product Information

Construct:	bFGF(134-288)
Species:	Human
Lyophilized From:	0.2 μ m filtered solution containing 5 mM Tris, pH 7.5, 150 mM NaCl
Expression System:	<i>E. coli</i>
Format:	Lyophilized powder
Stability:	Stable for at least 2 years from date of receipt at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
Genbank Accession:	NM_002006
MW:	17 kDa
Purity:	>98% by SDS-PAGE and HPLC analysis
Endotoxin Level:	< 1EU per ug of bFGF as determined by the LAL method
Biological Activity:	Measured by its ability to stimulate proliferation of NIH3T3 cells. The ED50 of this effect is between 0.1 - 1 ng/ml.
Reconstitution:	Reconstitute at 0.1-1.0 mg/ml in distilled water. This solution can then be diluted into other buffers. To maximize product collection from the vial surface, vortex briefly and then spin down to recollect the liquid.
Applications:	Useful for cell culture and for the study of signaling pathways.