

Recombinant Human FGF2

Recombinant human protein expressed in E. coli cells

Catalog # F820-40N

Lot # G3211-34

Product Description

Recombinant Human FGF2 was expressed in E. coli cells. The protein accession number is [P09038](#)

Gene Aliases

bFGF, Fibroblast growth factor 2, FGF-2, Heparin-binding growth factor 2, HBGF-2
Recombinant Human Basic Fibroblast Growth Factor (FGF2)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered solution in PBS

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in sterile distilled water to a concentration not less than 0.1 mg/mL is recommended. Please note, filter sterilization is a must following reconstitution. This solution can then be diluted into other buffers.

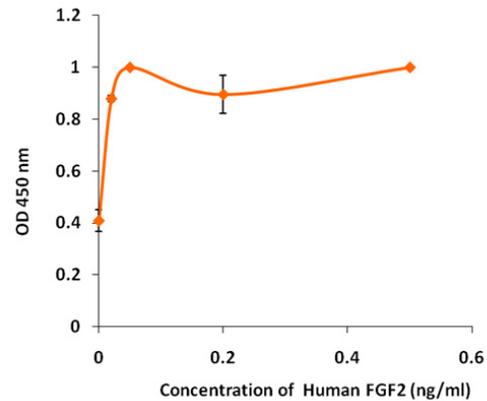
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

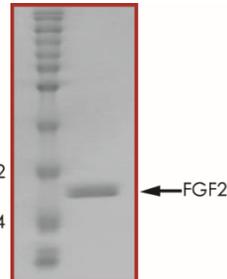
Basic Fibroblast Growth Factor is found in almost all tissues of mesodermal and neuroectodermal origin and also in tumors derived from these tissues. Endothelial cells produce large amounts of this factor. Some bFGF is associated with the extracellular matrix of the subendothelial cells. Many cells express bFGF only transiently and store it in a biologically inactive form. The mechanism by which the factor is released by the cells is not known. It is released after tissue injuries and during inflammatory processes. FGF receptors are encoded by a gene family consisting of at least four receptor tyrosine kinases that transduce signals important in a variety of developmental and physiological processes related to cell growth and differentiation. bFGF stimulates the growth of fibroblasts, myoblasts, osteoblasts, neuronal cells, endothelial cells, keratinocytes, chondrocytes, and many other cell types.

Specific Activity



The bioactivity was determined in a NIH/3T3 cell proliferation assay. The ED50 was in the range of 0.05 - 0.5 ng/ml.

Purity



Recombinant Human FGF2 resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.0 kDa**

Recombinant Human FGF2

Recombinant human protein expressed in E. coli cells

Catalog Number	F820-40N
Specific Lot Number	G3211-34
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.