

Recombinant Human KGF (FGF7)

Recombinant human protein expressed in E. coli cells

Catalog # K818-40N

Lot # G3211-84

Product Description

Recombinant Human KGF (FGF7) was expressed in E. coli cells. The protein accession number is [P21781](#)

Gene Aliases

KGF, Fibroblast Growth Factor-7, HBGF-7, Heparin-binding growth factor 7, Recombinant Human Keratinocyte Growth Factor (FGF7)

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 (pH 7.0)

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

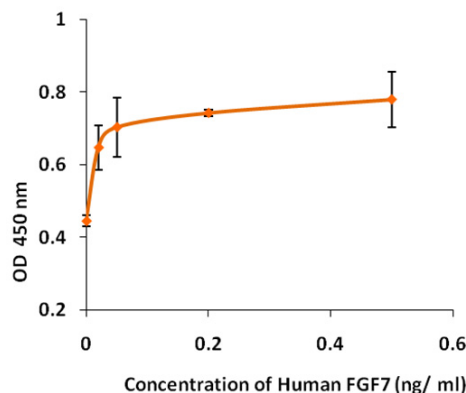
Storage and Stability

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

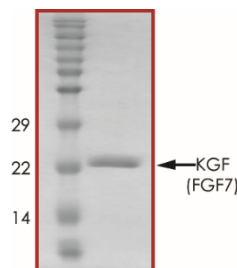
KGF is a member of the family of fibroblast growth factors, it is secreted in large amounts by fibroblast-like stromal cells in epithelial tissues. The KGF receptor encodes a tyrosine kinase and is a member of the family of receptors binding aFGF and bFGF. Binding of KGF to its receptor is competed by aFGF and bFGF. Although KGF is a potent specific mitogen for many epithelial cells it is not so for fibroblasts and endothelial cells. KGF is thought to play an important role in the paracrine growth control of normal epithelial cells. KGF stimulates the proliferation of primary and secondary human keratinocytes to the same extent as EGF.

Specific Activity



The ED50 as determined by the dose-dependent proliferation of 3T3 cells was found to be less than 0.5 ng/mL.

Purity



Recombinant Human KGF (FGF7) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **19.0 kDa**

Recombinant Human KGF (FGF7)

Recombinant human protein expressed in E. coli cells

Catalog Number	K818-40N
Specific Lot Number	G3211-84
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.