

Catalog #	Aliquot Size
F840-40N-30	30 µg
F840-40N-150	150 µg
F840-40N-1000	1 mg

Recombinant Human FGF20

Recombinant human protein expressed in E. coli cells

Catalog # F840-40N

Lot # G3211-35

Product Description

Recombinant Human FGF20 was expressed in E. coli cells. The protein accession number is [Q9NP95](#)

Gene Aliases

FGF-20, Recombinant Human Fibroblast Growth Factor 20 (FGF20)

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 10mM MOPS, 10mM ammonium sulphate, 0.2mM DTT and 0.2mM EDTA (pH 7.0).

Reconstitution Protocol

Reconstitute at 10 µg/mL in sterile PBS containing at least 0.1% human or bovine serum albumin

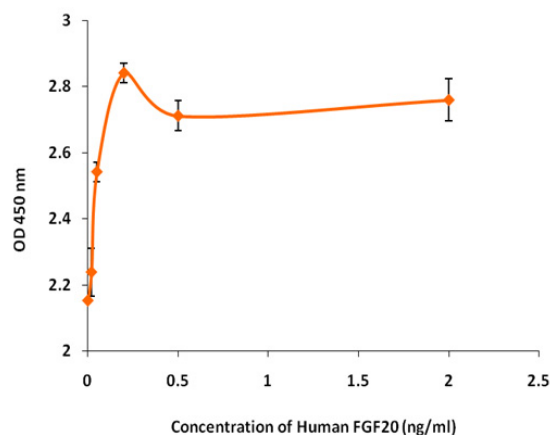
Storage and Stability

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

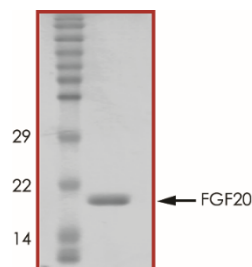
FGF-20 is a heparin binding protein of the FGF family which works to control cell division, differentiation, and function. These proteins are especially prominent during prenatal development and postnatal growth. FGF-20 shares a high degree of homology with FGF-9 and FGF-16. The group of receptors which FGF-20 binds to comprises of FGFR1c, FGFR2c, FGFR3b, FGFR3c and FGFR4. It is shown that FGF-20 is found in a variety of cells, including but limited to, dopaminergic neurons, fibroblasts, keratinocytes and breast epithelium.

Specific Activity



Measured in a cell proliferation assay using NR6R-3T3 mouse fibroblast cells

Purity



Recombinant Human FGF20 resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **23.5 kDa**

Recombinant Human FGF20

Recombinant human protein expressed in E. coli cells

Catalog Number	F840-40N
Specific Lot Number	G3211-35
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.