

FGF-basic 154 (human), (recombinant)

Fibroblast growth factors (FGFs) are a 22 member family of proteins known to be involved in angiogenesis, wound healing and embryonic development. As a family, they bind to heparin and signal through four receptor tyrosine kinases called, FGFR1, 2, 3 and 4. Although the mechanism remains unclear, FGF-basic 154, also known as FGF-2, is a critical component in keeping embryonic stem cells undifferentiated in cell culture systems.

Ordering Information [Order Online »](#)

ALX-201-823-0050	50µg
------------------	------

Manuals, SDS & CofA [View Online »](#)

- Carrier-free

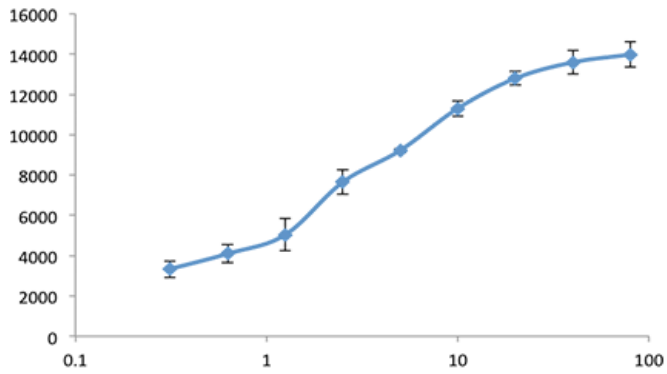


Figure 1: Serial dilutions of Human FGF Basic 154, starting at 80 ng/mL, were added to NIH 3T3 cells. Cell proliferation was measured after 48 hours and the linear portion of the curve can be used to calculate the ED50.

Handling & Storage

Use/Stability	It is recommended that a carrier protein (0.1% HSA or BSA) is added for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	FGF2, HBGF-2, Prostatropin
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the dose-dependent proliferation of mouse BALB/c 3T3 cells which is typically observed at concentrations <1ng/ml.
Endotoxin Content	≤1 EU/μg protein measured by kinetic LAL analysis.
Formulation	Lyophilized from 10mM sodium phosphate, 75mM NaCl, pH 7.5. Sterile filtered.
MW	~17.2kDa
Purity	≥95% (Reducing and Non-reducing SDS-PAGE)
Reconstitution	When reconstituting the product, gently pipet and wash down the sides of the vial to ensure full recovery of the protein into solution. It is recommended to reconstitute the lyophilized product with sterile water at a concentration of 0.1 mg/ml, which can be further diluted into other aqueous solutions.
Source	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 154 amino acids.
UniProt ID	P09038 + 7 aa of propeptide

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)