

# FGF-basic 147 (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 147 (variant of FGF basic 154), also called FGF-2, is a critical component in keeping embryonic stem cells undifferentiated in cell culture systems.

## Ordering Information

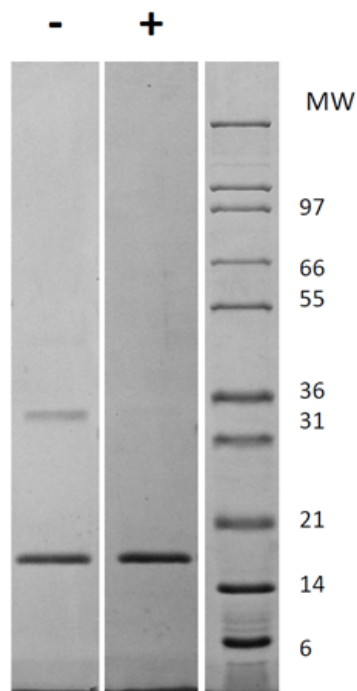
[Order Online »](#)

ALX-201-822-0010	10µg
ALX-201-822-0050	50µg
ALX-201-822-0100	100µg

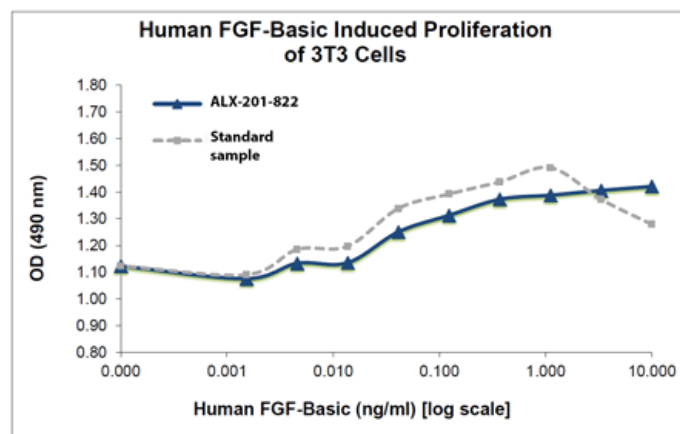
## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 2:** 1 $\mu$ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human FGF basic has a predicted MW of 16.5 kDa.



**Figure 1: Human FGF-basic 147 Bioactivity Data.** 3T3 cells were cultured with 0 to 10ng/ml human FGF basic. Cell proliferation was measured after 44 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of human FGF basic was 0.042-0.063ng/ml. The typical expected range is less than 1ng/ml.

# Handling & Storage

Use/Stability	It is recommended that working aliquots to be diluted in a 0.1% BSA solution for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -80°C and avoid repeat freeze thaws.
Long Term Storage	-80°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	≤5 ng/mL. The ED50 determined by NR6R-3T3 cell proliferation assay
Endotoxin Content	≤0.1 EU/μg using USP /EP 2.6.14
Formulation	Lyophilized from 10mM sodium phosphate, 75mM sodium chloride, pH 7.5. Sterile filtered.
MW	~16.5kDa
Purity	≥95% using Coomassie Staining
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 147 amino acids.
UniProt ID	P09038



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)