

# FGF basic (mouse), (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, or FGF-1, is a critical component in keeping human embryonic stem cells undifferentiated in cell culture systems.

## Ordering Information

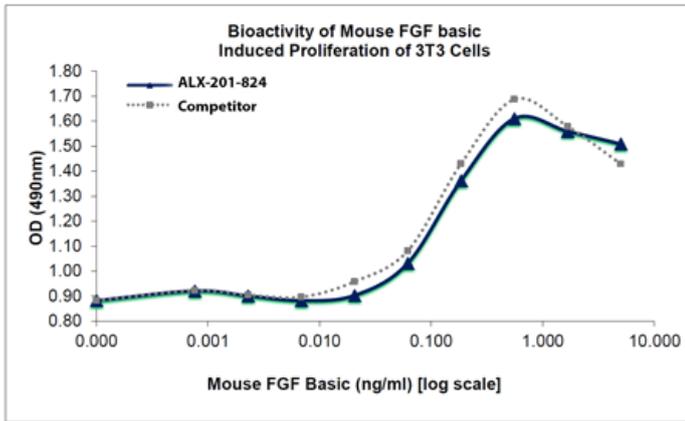
[Order Online »](#)

ALX-201-824-0050	50µg
------------------	------

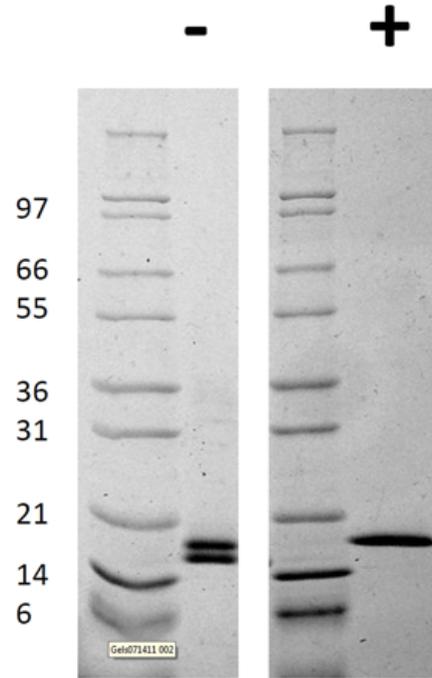
## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 1: Mouse FGF basic Bioactivity Data.** Serial dilutions of mouse FGF Basic, starting at 5ng/ml, were added to NIH 3T3 cells. Cell proliferation was measured after 41 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of mouse FGF Basic was 0.1-0.16ng/ml. This value is comparable with the typical expected range of < 1ng/ml.



**Figure 2:** 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Mouse FGF-basic has a predicted MW of 17.2 kDa.

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	FGF2, HBGF-2, Prostatropin
<b>Appearance</b>	White lyophilized (freeze-dried) powder.
<b>Biological Activity</b>	The activity is determined by the dose-dependent proliferation of BALB/3T3 cells which is typically observed at concentrations <1ng/ml.
<b>Endotoxin Content</b>	≤1 EU/μg protein measured by kinetic LAL analysis.
<b>Formulation</b>	Lyophilized from a 0.2 μm filtered solution containing 10 mM Sodium Phosphate, 50 mM Sodium Chloride, pH 7.5.
<b>MW</b>	~16.3kDa
<b>Purity</b>	≥95% (Reducing and Non-reducing SDS-PAGE)
<b>Reconstitution</b>	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.
<b>Source</b>	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 145 amino acids.
<b>UniProt ID</b>	p15655

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)