

# FGF-5 (human), (recombinant)

FGF-5 functions as a regulatory factor during hair elongation and skeletal muscle development. Fibroblast growth factor 5 (FGF-5) is a secreted heparin-binding growth factor that binds to FGF receptors 1 and 2 (FGFR1 and FGFR2).

## Ordering Information

[Order Online »](#)

ENZ-PRT231-0050	50µg
ENZ-PRT231-0100	100µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Fibroblast growth factor 5, Heparin-binding growth factor 5, HBGF-5, Smag-82
Biological Activity	3T3 proliferation with 1 µg heparin; ED50 ≤ 10 ng/mL ( $\geq 1.0 \times 10^5$ units/mg).
Endotoxin Content	≤1 EU/µg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium chloride, pH 7.5.
MW	~27.7 kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 252 amino acids.
UniProt ID	P12034



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)