

# FGF-acidic (human), (recombinant)

Acidic fibroblast growth factor (FGF-acidic), is a potent inducer of DNA synthesis, cell proliferation, and has chemotactic activities. FGF-acidic regulates cardiogenesis through protein kinase C signaling. FGF-acidic also functions as an insulin sensitizer and mediates adipose tissue remodeling.

## Ordering Information

[Order Online »](#)

ENZ-PRT235-0050	50µg
ENZ-PRT235-0100	100µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Fibroblast growth factor 1, FGF-1, HBGF-1,  $\beta$ -endothelial growth factor, ECCF- $\beta$

**Biological Activity** 3T3 proliferation; ED50  $\leq 2$  ng/mL ( $\geq 5.0 \times 10^5$  units/mg).

**Endotoxin Content**  $\leq 1$  EU/ $\mu$ g protein (LAL test)

**Formulation** Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium phosphate, 150 mM sodium sulfate, pH 7.5.

**MW** ~16.0 kDa

**Purity**  $\geq 95\%$  (SDS-PAGE)

**Reconstitution** Reconstitute in sterile water at 0.1 mg/mL.

**Source** Produced in *E. coli*. Contains 141 amino acids.

**UniProt ID** P05230



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-  
fr@enzolifesciences.com](mailto:info-fr@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)