

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.

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, β -endothelial growth factor, ECCF- β
Biological Activity	3T3 proliferation; ED50 \leq 2 ng/mL (\geq 5.0 x 10 ⁵ units/mg).
Endotoxin Content	\leq 1 EU/ μ 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 <i>E. coli</i> . Contains 141 amino acids.
UniProt ID	P05230

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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