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 NameFibroblast 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 10⁵ units/mg).

Endotoxin Content ≤1 EU/μg 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 ≥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

Last modified: May 29, 2024

