# FGF-6 (human), (recombinant)

Fibroblast growth factor 6 (FGF-6) is an important factor driving muscle differentiation and regeneration. FGF-6 is a heparin-binding growth factor that is expressed in epithelial and mesenchymal lineages. FGF-6 binds and signals through the FGF receptors FGFR1, FGFR2, and FGFR4. FGF-6 functions as a mitogen for vascular endothelial cells and fibroblasts.

## **Ordering Information**

Order Online »

ENZ-PRT232-0025	25µg
ENZ-PRT232-0100	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 6, Heparin secretory-transforming

protein 2, HST-2, HSTF-2, HBGF-6

Biological Activity 3T3 proliferation with 1 µg heparin; ED50 ≤ 1 ng/mL (≥ 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 phosphate and 50 mM sodium

chloride, pH 7.5.

**MW** ~18.9 kDa

Purity ≥95% (SDS-PAGE)

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

**Source** Produced in *E. coli*. Contains 169 amino acids.

UniProt ID P10767

Last modified: May 29, 2024

