

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.

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 x 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 <i>E. coli</i> . Contains 169 amino acids.
UniProt ID	P10767

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