

Data Sheet

CellGenix® Recombinant Human Fibroblast Growth Factor-2 (rh FGF-2)

Preclinical Grade - Order No.: 1421-010 (10 µg)*, 1421-050 (50 µg)

Product Characteristics

Source	<i>E. coli</i>
Description	Human FGF-2, accession # P09038, Ala135-Ser288 N-terminal Met Molecular mass 17.3 kDa
Formulation	Lyophilized from a 0.2 µm-filtered solution containing 25 mM sodium phosphate, 400 mM sodium chloride, 10 mM glutathione (red.), and 5 % mannitol, pH 6.5
Intended use	For preclinical ex vivo use. Not intended for therapeutic use.

Quality Parameters

Activity	≥ 0.9 x 10 ⁶ IU/mg calibrated against NIBSC #90/712 Measured in a cell proliferation assay using an FGF-2-dependent cell line, FBHE
Purity	≥ 95 %, as determined by SDS-PAGE (under reducing conditions, visualized by Coomassie staining) and RP-HPLC
Endotoxin	< 1000 EU/mg, as determined by LAL gel clot test
Sterility	Sterility test of the vial product
Mass per vial	1421-010: 10 µg, 1421-050: 50 µg
Animal-derived component-free	ADCF Level 1: The final product contains neither animal- nor human-derived materials.

Shipment & Storage

Transport	Ambient temperature. Please refer to Technote (www.cellgenix.com)
Shelf life	3 years from date of shipment
Storage & Stability	Store lyophilized cytokine at -20 °C to -80 °C. <ul style="list-style-type: none">• 4 months at -20 °C to -80 °C under sterile conditions after reconstitution. Store in aliquots in polypropylene cryogenic vials. Avoid repeated freeze/thaw cycles.

Handling Instructions

Reconstitution	Recommended in sterile water to a final concentration of 100 µg/ml for 10 µg vials or 250 µg/ml for 50 µg vials.
Dilution	Recommended in CellGenix® serum-free media. For dilution with protein free medium, a carrier protein (0.1-1 % albumin or 1-10 % appropriate serum) has to be included. Failure to dilute product according to these instructions may result in loss of activity.

Quality Statement

Final manufacturing steps and QC are performed in a GMP facility. No animal- or human-derived components are present in the final product (ADCF Level 1).