

## Data Sheet

# CellGenix® Recombinant Human Epidermal Growth Factor (rh EGF) Preclinical Grade – Order No.: 1416-050 (50 µg)

## Product Characteristics

Source	<i>E. coli</i>
Description	Human EGF, accession # P01133, Asn971-Arg1023 N-terminal Met and C-terminal 6xHis-tag Molecular mass 7.2 kDa
Formulation	Lyophilized from a 0.2 µm-filtered solution containing 25 mM sodium acetate, pH 5.0
Intended use	For preclinical ex vivo use. Not intended for therapeutic use.

## Quality Parameters

Activity	≥ 0.5 x 10 <sup>6</sup> IU/mg calibrated against NIBSC #91/530 Measured in a cell proliferation assay using an EGF-dependent cell line, Balb/3T3
Purity	≥ 95 %, as determined by RP-HPLC
Endotoxin	< 1000 EU/mg, as determined by LAL gel clot test
Sterility	Sterility test of the vial product
Mass per vial	50 µg
Animal-derived component-free	<b>ADCF Level 2:</b> The final product contains neither animal- nor human-derived materials. ADCF Level 2 cytokines are produced in our dedicated animal-free facility. No animal-derived components are used throughout the complete production process. All ADCF Level 2 cytokines are produced in <i>E. coli</i> .

## Shipment & Storage

Transport	Ambient temperature. Please refer to Technote ( <a href="http://www.cellgenix.com">www.cellgenix.com</a> )
Shelf life	3 years from date of shipment
Storage & Stability	Store lyophilized cytokine at -20 °C to -80 °C. Store a 250 µg/ml cytokine solution: <ul style="list-style-type: none"><li>• 4 weeks at 2 °C to 8 °C under sterile conditions after reconstitution. Store in the original container.</li><li>• 4 months at -20 °C to -80 °C under sterile conditions after reconstitution. Store in aliquots in polypropylene cryogenic vials.</li></ul> Avoid repeated freeze/thaw cycles.

## Handling Instructions

Reconstitution	Recommended in 0.2 % acetic acid to a final concentration of 250 µg/ml.
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 and no animal- or human-derived materials were used in production (ADCF Level 2).