

# CellGenix® GMP Recombinant Human Epidermal Growth Factor (rh EGF)



## Product Information

CellGenix® Recombinant Human EGF reliably promotes the expansion of mesenchymal stem cells (MSCs). In addition, it promotes the differentiation of embryonic stem cells (ESCs) and induced pluripotent stem cells (iPSCs)-derived hepatoblasts. CellGenix® rh EGF is produced in our dedicated animal-free facility ensuring maximum safety for optimal use in ATMP manufacturing. 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).

## Features and Benefits

- Seamless transition from early development to clinical stages with consistent product quality & performance
- Exceptional batch-to-batch consistency to ensure optimal cell expansion and differentiation
- Manufactured in compliance with applicable GMP guidelines and in accordance with USP Chapter <1043>, Ph. Eur. General Chapter 5.2.12 and ISO 20399:2022
- We offer expert regulatory support in all phases from development to post-approval to assist you in safely bringing your product to the market. Drug Master File filed FDA CBER (USA)

## Application

Epidermal Growth Factor (EGF) is a growth factor that promotes cell growth, proliferation and differentiation. It is found in a variety of human body fluids.

EGF is used in the cell and gene therapy space for the expansion of bone marrow and adipose tissue derived MSCs, and the generation of stem cell derived hepatocytes.

## Product Characteristics

<b>Source</b>	<i>E. coli</i>
<b>Description</b>	Human EGF, accession # P01133, Asn971-Arg1023 N-terminal Met and C-terminal 6xHis-tag Molecular mass 7.2 kDa
<b>Formulation</b>	Lyophilized from a 0.2 µm-filtered solution containing 25 mM sodium acetate, pH 5.0
<b>Intended use</b>	For further manufacturing use.

## Quality Parameters

<b>Identity</b>	Confirmed by DNA-sequencing of the expression plasmid in end-of-production cells and N-terminal sequencing of the final product
<b>Activity</b>	0.5 – 2.0 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
<b>Purity</b>	≥ 97%, as determined by RP-HPLC
<b>Product-related proteins</b>	≤ 5% as determined by RP-HPLC
<b>Host-cell DNA</b>	≤ 200 ng/mg, as determined with a fluorimetric assay
<b>Host-cell protein</b>	≤ 1.0 µg/mg, as determined by ELISA
<b>Endotoxin</b>	≤ 50 EU/mg, as determined by LAL gel clot test according to Ph. Eur. and USP
<b>Sterility</b>	Sterility test according to Ph. Eur. and USP of the vial product
<b>Mass per vial</b>	<b>50 µg:</b> 43 - 57 µg, <b>1000 µg:</b> 900 – 1100 µg as determined by spectrophotometrical measurement
<b>Animal-derived component-free</b>	<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 and Storage

<b>Transport</b>	Ambient temperature. Please refer to Technical Note "Shipment of CellGenix® Preclinical and GMP Cytokines at Ambient Temperatures".
<b>Shelf Life</b>	Minimum 6 months from date of shipping
<b>Storage and Stability</b>	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

<b>Reconstitution</b>	Recommended in 0.2% acetic acid to a final concentration of 250 µg/mL for 50 µg vials or 500 µg/mL for 1000 µg vials.
<b>Dilution</b>	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.

## Packaging

CellGenix® cytokines are provided in glass vials, closed with vacuum rubber stoppers and sealed with aluminum tear off caps. The following material is used:

### Glass vials

For 50 µg vials: Glass vials (2 mL; colorless; 35.00 x 13.75 mm) with DIN Crimp Neck N13-2 made from borosilicate glass hydrolytic type I (in compliance with Ph. Eur. 3.2.1 and USP <660> glass containers for pharmaceutical use).

For 1 mg vials: Glass vials (6 mL; colorless; 40.00 x 22.00 mm) with DIN Crimp Neck N20 made from borosilicate glass hydrolytic type I (in compliance with Ph. Eur. 3.2.1 and USP <660> glass containers for pharmaceutical use).

### Vacuum rubber stoppers, Type I butyl rubber

The formulation is 4023/50/grey. This corresponds to bromobutyl rubber with a hardness of 50 (hardness measured in shore A). This is compliant with Ph. Eur. 3.2.9 Type 1 and with the physicochemical tests as described in USP General Chapter <381> "Elastomeric Closures for Injections".

### Aluminum tear off caps

Aluminum tear off caps (13 mm; gold) are produced in accordance to valid quality criteria for metal caps.

The container closure has been validated after a storage period of up to 5 years at -80 °C by verification of sterility. In addition, the container closure has been demonstrated according to USP <671>.

## Ordering Information

Product Description	Size & Package	Storage	Cat. No.
CellGenix® GMP EGF	50 µg	-20 °C to -80 °C	1016-050
CellGenix® GMP EGF	1 mg	-20 °C to -80 °C	1016-1000

## Sartorius is Your Reliable Supply Partner

High-quality raw materials are essential to ensure safety, efficacy and batch-to-batch consistency. We propose premium-grade raw materials suitable from preclinical development to the manufacturing of the therapy. Our GMP grade products allow for the safe use in clinical trials and commercial manufacturing.

Our GMP cytokines include documented evidence of lot specific sterility, activity, and shelf-life. Our experts will help simplify your raw material qualification and validation efforts. We provide customized solutions to your enquiries, as well as quality control services to ensure the quality of our products. Our regulatory expertise guarantees a suited service to your regulatory procedures, ensuring an extensive support every step of the process.

## Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

Sartorius CellGenix GmbH  
Am Flughafen 16  
79108 Freiburg  
Phone +49 761 88889 0  
Fax + 49 761 88889 830  
info-freiburg@sartorius.com

## USA

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

 **For more information, visit**  
[sartorius.com](https://www.sartorius.com)