

Data Sheet

CellGenix® GMP Recombinant Human Interleukin-3 (rh IL-3)

Order No.: 1002-050 (50 µg)

Product Characteristics

| | |
|--------------|--|
| Source | <i>E. coli</i> |
| Description | Human Interleukin-3, accession # P08700, Ala20-Phe152 N-terminal Met Molecular mass 15.2 kDa |
| Formulation | Lyophilized from a 0.2 µm-filtered solution containing 10mM sodium phosphate and 50 mM sodium chloride, pH 7.2 |
| Intended use | For further manufacturing use. |

Quality Parameters

| | |
|-------------------------------|---|
| Identity | Confirmed by DNA-sequencing of the expression plasmid in end-of-production cells and N-terminal sequencing of the final product |
| Activity | 0.9 – 3.6 x 10 ⁶ IU /mg calibrated against NIBSC #91/510 Measured in a cell proliferation assay using an IL-3-dependent cell line, TF-1 |
| Purity | ≥ 97 %, as determined by SDS-PAGE (under reducing and non-reducing conditions, visualized by Coomassie staining) and RP-HPLC |
| Product-related proteins | ≤ 5 % oligomers as determined by SDS-PAGE (under non-reducing conditions, visualized by Coomassie staining) and ≤ 20 % oxidized IL-3 species as determined by RP-HPLC |
| Host-cell DNA | ≤ 200 ng/mg, as determined with a fluorimetric assay |
| Host-cell protein | ≤ 1.0 µg/mg, as determined by ELISA |
| Endotoxin | ≤ 50 EU/mg, as determined by LAL gel clot test according to Ph. Eur. and USP |
| Sterility | Sterility test according to Ph. Eur. and USP of the vialled product |
| Mass per vial | 43 – 57 µg, as determined by spectrophotometrical measurement |
| Animal-derived component-free | ADCF Level 2: 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 (www.cellgenix.com) |
| Shelf life | Minimum 6 months from date of shipping |
| Storage & Stability | Store lyophilized cytokine at -20 °C to -80 °C. <ul style="list-style-type: none">• Store a 250 µg/ml reconstituted cytokine solution for 4 weeks at 2 °C to 8 °C under sterile conditions after reconstitution. Store in the original container.• Store a 100 µg/ml reconstituted cytokine solution for 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 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

This product is manufactured, tested and released in compliance with the relevant GMP-guidelines. 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). USP chapter <1043> has been considered in the design of this product. This product is compliant with Ph. Eur. general chapter 5.2.12 and ISO 20399:2022.