

## Data Sheet

### CellGenix® Recombinant Human Interleukin-3 (rh IL-3) Preclinical Grade – Order No.: 1402-010 (10 µg)\*, 1402-050 (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 10 mM sodium phosphate and 50 mM sodium chloride, pH 7.2.
Intended use	For preclinical ex vivo use. Not intended for therapeutic use.

#### Quality Parameters

Activity	≥ 0.9 x 10 <sup>6</sup> IU/mg calibrated against NIBSC #91/510 Measured in a cell proliferation assay using an IL-3-dependent cell line, TF-1
Purity	≥ 95 %, as determined by SDS-PAGE (under reducing and non-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	1402-010: 10 µg, 1402-050: 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

<b>Transport</b>	Ambient temperature. Please refer to Technote ( <a href="http://www.cellgenix.com">www.cellgenix.com</a> )
<b>Shelf life</b>	3 years from date of shipment
<b>Storage &amp; Stability</b>	<p>Store lyophilized cytokine at -20 °C to -80 °C.</p> <ul style="list-style-type: none"><li>• 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.</li><li>• 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.</li></ul> <p>Avoid repeated freeze/thaw cycles.</p>

## Handling Instructions

<b>Reconstitution</b>	Recommended in sterile water to a final concentration of 100 µg/ml (for 10 µg vials) or 250 µg/ml (for 50 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.

## 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).