

## Data Sheet

### CellGenix® Recombinant Human Interleukin-2 (rh IL-2) Preclinical Grade – Order No.: 1420-010 (10 µg)\*, 1420-050 (50 µg)

#### Product Characteristics

Source	<i>E. coli</i>
Description	Human Interleukin-2 with a substitution of cysteine-125 by serine, accession # P60568, Pro22–Thr153 N-terminal Met and C-terminal 6xHis-tag Molecular mass 16.3 kDa With the exception of the added C-terminal 6xHis-tag, the amino acid sequence corresponds to that of Aldesleukin
Formulation	Lyophilized from a 0.2 µm-filtered solution containing 25 mM sodium acetate and 50 mM sodium chloride, pH 5.0
Intended use	For preclinical <i>ex vivo</i> use. Not intended for therapeutic use.

#### Quality Parameters

Activity	≥ 8 x 10 <sup>6</sup> IU/mg calibrated against an internal reference standard* Measured in a cell proliferation assay using an IL-2-dependent cell line, CTLL-2. <i>*The internal reference standard was calibrated against Proleukin®. Proleukin® is a registered trademark of Novartis AG</i>
Purity	≥ 95 %, as determined by SDS-PAGE (under reducing and non-reducing conditions, visualized by Coomassie staining) and RP-HPLC
Endotoxin	< 25 EU/mg, as determined by LAL gel clot test
Sterility	Sterility test of the vial product
Mass per vial	1420-010: 10 µg, 1420-050: 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>	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 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).