

Data Sheet

CellGenix® Recombinant Human Interleukin-7 (rh IL-7) Preclinical Grade – Order No.: 1410-010 (10 µg)*, 1410-050 (50 µg)

Product Characteristics

Source	<i>E. coli</i>
Description	Human Interleukin-7, accession # P13232, Asp26-His177 N-terminal Met and C-terminal 6xHis-tag Molecular mass 18.3 kDa
Formulation	Lyophilized from a 0.2 µm-filtered solution containing 25 mM sodium phosphate, 150 mM sodium chloride, and 0.1% polysorbate 20, pH 7.5.
Intended use	For preclinical ex vivo use. Not intended for therapeutic use.

Quality Parameters

Activity	≥ 50 x 10 ⁶ IU/mg calibrated against NIBSC #90/530 Measured in a cell proliferation assay using an IL-7-dependent cell line, IxN/2b
Purity	≥ 95 %, as determined by SDS-PAGE (under reducing and non-reducing conditions, visualized by Coomassie staining)
Endotoxin	< 25 EU/mg, as determined by LAL gel clot test
Sterility	Sterility test of the vial product
Mass per vial	1410-010: 10 µg, 1410-050: 50 µg
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	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">• 4 weeks at 2 °C to 8 °C under sterile conditions after reconstitution. Store in the original container.• 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 100 µg/ml (for 10 µg vials) or 250 µg/ml (for 50 µg vials).
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).