

## Data Sheet

# CellGenix® Recombinant Human Interferon-gamma (rh IFN- $\gamma$ )

Preclinical Grade – Order No.: 1425-050 (50  $\mu$ g)

### Product Characteristics

Source	<i>E. coli</i>
Description	Human Interferon-gamma, accession # P01579, Gln24-Arg162 N-terminal Met Molecular mass 16.5 kDa
Formulation	Lyophilized from a 0.2 $\mu$ m-filtered solution containing 20 mM succinate, pH 5.0.
Intended use	For preclinical <i>ex vivo</i> use. Not intended for therapeutic use.

### Quality Parameters

Activity	$\geq 10 \times 10^6$ IU/mg calibrated against Imukin® Measured in a cell cytotoxicity assay using a IFN- $\gamma$ -sensitive cell line, HT-29
Purity	$\geq 95$ %, as determined by SDS-PAGE (under reducing and non-reducing conditions, visualized by Coomassie staining)
Endotoxin	$< 1000$ EU/mg, as determined by LAL gel clot test
Sterility	Sterility test of the vial product
Mass per vial	50 $\mu$ g
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

Transport	Ambient temperature. Please refer to Technote ( <a href="http://www.cellgenix.com">www.cellgenix.com</a> )
Shelf life	3 years from date of shipment
Storage & Stability	Store lyophilized cytokine at -20°C to -80°C. 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

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).