

## Data Sheet

CellGenix® GMP Dendritic Cell Medium (DC), Serum-free, Xeno-free  
 Order No.: 20801-0500 (500 ml bottle), 20901-0500 (500 ml bag)

### Product Characteristics

<b>Main components</b>	Salts, sugars, amino acids, vitamins, buffers Contains phenol red and L-glutamine
<b>Human proteins</b>	<b>Transferrin:</b> human plasma-derived. <b>Albumin:</b> the plasma used in the manufacture of HSA complies with the requirements of the Ph. Eur. Monograph "Human Plasma for Fractionation" (0853) and European regulations. The HSA is manufactured, tested and released according to the current GMP requirements and the current Ph. Eur. The HSA manufacturer complies with directive 2002/98/EC setting standards of quality and safety for the collection, testing, processing, storage and distribution of human blood and blood components. <b>Insulin:</b> human recombinant, yeast-derived, complies with Ph. Eur. and USP.
<b>Application</b>	Generation of dendritic cells
<b>Intended use</b>	For further manufacturing use.

### Quality Parameters

<b>Appearance</b>	Clear, red liquid
<b>pH</b>	7.2 - 7.5, determined according to Ph. Eur.
<b>Osmolality</b>	260 - 320 mOsm/kg H <sub>2</sub> O, determined according to Ph. Eur.
<b>Bioassay</b>	Generation of human dendritic cells from CD14 <sup>+</sup> monocytes
<b>Endotoxin</b>	≤ 1 EU/ml, determined according to Ph. Eur.
<b>Mycoplasma</b>	Not detectable, determined according to Ph. Eur.
<b>Sterility</b>	Sterility test of the final product, determined according to Ph. Eur.

### 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 production. Minimum 6 months from date of shipping.
<b>Storage</b>	Store at +2 °C to +8 °C. Light protection is recommended.

## Quality Statement

This product is manufactured, tested and released in compliance with the relevant GMP-guidelines. USP chapter <1043> has been considered in the design of this product. This product is compliant with ISO 20399:2022. The formulation does not contain animal-derived components (xeno-free).

Human proteins have been collected from healthy donors at the time of collection, and samples of their donations were tested individually and found negative for viral diseases by approved methods (HIV1/HIV2, HBV, HCV, Parvovirus B19).