

SECTION 1: Identification of the substance/mixture and of the company/undertaking**1.1. Product identifier**

| | |
|--------------|--|
| Product form | : Mixture |
| Trade name | : CellGenix® GMP Dendritic Cell Medium (DC), Serum-free, w/o Phenol Red |
| Product code | : 20805 (bottle), 20905 (bag) |

1.2. Relevant identified uses of the substance or mixture and uses advised against**1.2.1. Relevant identified uses**

Use of the substance/mixture : Cell Culture component. For further manufacturing use.

1.2.2. Uses advised against

Restrictions on use : Not intended for human in vivo application.

1.3. Details of the supplier of the safety data sheet**Supplier**

Sartorius CellGenix GmbH
Am Flughafen 16
79108 Freiburg
T +49 761 888 89-0
F +49 761 888 89-830

E-mail address of competent person responsible for the SDS

info@ubsplus.de

1.4. Emergency telephone number

Emergency number : +49 761 19240
(VIZ Freiburg, 24 h, German & English)

SECTION 2: Hazards identification**2.1. Classification of the substance or mixture**

Classification according to Regulation (EC) No. 1272/2008 [CLP] and 29 CFR § 1910.1200, Appendix A and B

Not classified

2.2. Label elements

Labelling according to Regulation (EC) No. 1272/2008 [CLP] and 29 CFR § 1910.1200, Appendix C

No labelling applicable

2.3. Other hazards

This substance/mixture does not meet the PBT criteria of REACH regulation, annex XIII

This substance/mixture does not meet the vPvB criteria of REACH regulation, annex XIII

SECTION 3: Composition/information on ingredients**3.1. Substances**

Not applicable

3.2. Mixtures

This mixture does not contain any substances to be mentioned according to the criteria of section 3.2 of REACH annex II

SECTION 4: First aid measures**4.1. Description of first aid measures**

| | |
|-------------------------------------|---|
| First-aid measures general | : Move the affected person away from the contaminated area. Fresh air, rest. Prevent cooling by covering the victim (no warming up). If unconscious place in recovery position and seek medical advice. Do not give an unconscious person anything to drink. Remove soiled clothing promptly. |
| First-aid measures after inhalation | : Remove person to fresh air and keep comfortable for breathing. Artificial respiration if indicated. If symptoms persist, call a physician. |

| | |
|---------------------------------------|---|
| First-aid measures after skin contact | : Immediately remove contaminated clothing or footwear. Rinse and then wash skin thoroughly with water and soap. If skin irritation or rash occurs: Get medical advice/attention. |
| First-aid measures after eye contact | : Wash with plenty of water (during 20 minutes minimum) with eyes wide open after taking off soft contact lenses and immediately take medical advice. |
| First-aid measures after ingestion | : Rinse mouth. Drink plenty of water. If you feel unwell, seek medical advice. |

4.2. Most important symptoms and effects, both acute and delayed

No additional information available

4.3. Indication of any immediate medical attention and special treatment needed

Treat symptomatically.

SECTION 5: Firefighting measures

5.1. Extinguishing media

| | |
|--------------------------------|--|
| Suitable extinguishing media | : water, carbon dioxide (CO ₂), powder and foam. Use extinguishing media appropriate for surrounding fire. |
| Unsuitable extinguishing media | : Not known. |

5.2. Special hazards arising from the substance or mixture

| | |
|--|---|
| Fire hazard | : The product is not flammable. |
| Explosion hazard | : Not known. |
| Hazardous decomposition products in case of fire | : Nitrogen oxides. Carbon monoxide. Carbon dioxide. |

5.3. Advice for firefighters

| | |
|--------------------------------|--|
| Firefighting instructions | : Do not allow run-off from fire fighting to enter drains or water courses. |
| Protection during firefighting | : Evacuate the danger area. Keep public away from danger area. Extra personal protection: complete protective clothing including self-contained breathing apparatus. |

SECTION 6: Accidental release measures

6.1. Personal precautions, protective equipment and emergency procedures

| | |
|------------------|--|
| General measures | : Do not breathe spray. Avoid contact with skin and eyes. Ventilate spillage area. In case of insufficient ventilation, wear suitable respiratory equipment. |
|------------------|--|

6.1.1. For non-emergency personnel

| | |
|----------------------|---|
| Protective equipment | : Avoid exposure. Wear personal protective equipment. |
|----------------------|---|

6.1.2. For emergency responders

| | |
|----------------------|---|
| Protective equipment | : Avoid exposure. Wear personal protective equipment. |
|----------------------|---|

6.2. Environmental precautions

Prevent entry to sewers and public waters.

6.3. Methods and material for containment and cleaning up

| | |
|-------------------------|---|
| For containment | : Dike and contain spill. Soak up spills with inert solids, such as clay or diatomaceous earth as soon as possible. |
| Methods for cleaning up | : Clean with water. |

6.4. Reference to other sections

No additional information available

SECTION 7: Handling and storage

7.1. Precautions for safe handling

| | |
|-----------------------------------|--|
| Additional hazards when processed | : Avoid formation of aerosols. Do not breathe vapour/aerosol. |
| Precautions for safe handling | : Avoid contact with skin, eyes and clothing. |
| Hygiene measures | : Do not eat, drink or smoke when using this product. Wash hands and other exposed areas with mild soap and water before eating, drinking or smoking and when leaving work. Wash contaminated clothing before reuse. |

7.2. Conditions for safe storage, including any incompatibilities

| | |
|---------------------|--|
| Storage conditions | : Keep container tightly closed. Protect from light. |
| Storage temperature | : 2 - 8 °C (Shipping at ambient temperature.) |

7.3. Specific end use(s)

Cell Culture component. For further manufacturing use. Not intended for human in vivo application.

SECTION 8: Exposure controls/personal protection

8.1. Control parameters

No additional information available

8.2. Exposure controls

| | |
|-----------------------------------|---|
| Appropriate engineering controls | : Ensure good ventilation of the work station. |
| Materials for protective clothing | : Wear proper protective equipment |
| Hand protection | : Chemically resistant protective gloves. Choosing the proper glove is a decision that depends not only on the type of material, but also on other quality features, which differ for each manufacturer. . Material : Latex (0,1 mm). Nitrile rubber (0,1 mm) |
| Eye protection | : Sealed safety goggles. (EN 166). |
| Respiratory protection | : No respiratory protection needed under normal use conditions |



Other information : Wash hands and other exposed areas with mild soap and water before eating, drinking or smoking and when leaving work.

SECTION 9: Physical and chemical properties

9.1. Information on basic physical and chemical properties

| | |
|--|------------------------------|
| Physical state | : Liquid |
| Colour | : Colourless |
| Odour | : odourless |
| Odour threshold | : Not applicable |
| pH | : 7.2 - 7.5 (Ph. Eur. 2.2.3) |
| pH solution | : Not determined |
| Relative evaporation rate (butylacetate=1) | : Not applicable |
| Melting point | : Not determined |
| Freezing point | : No data available |
| Boiling point | : Not determined |
| Flash point | : Not applicable |
| Auto-ignition temperature | : Not determined |
| Decomposition temperature | : Not determined |
| Flammability (solid, gas) | : Not applicable |
| Vapour pressure | : Not determined |
| Relative vapour density at 20 °C | : No data available |
| Relative density | : Not determined |
| Solubility | : Water: Soluble |
| Log Pow | : Not determined |
| Log Kow | : Not determined |
| Viscosity, kinematic | : Not determined |
| Viscosity, dynamic | : Not determined |
| Explosive properties | : Not known. |
| Oxidising properties | : Not known. |
| Explosive limits | : No data available |

9.2. Other information

No additional information available

SECTION 10: Stability and reactivity

10.1. Reactivity

No dangerous reactions known under normal conditions of use.

10.2. Chemical stability

Stable under normal conditions.

10.3. Possibility of hazardous reactions

No dangerous reactions known.

10.4. Conditions to avoid

No additional information available

10.5. Incompatible materials

Strong oxidizers.

10.6. Hazardous decomposition products

No additional information available

SECTION 11: Toxicological information

11.1. Information on toxicological effects

| | |
|-----------------------------------|---|
| Acute toxicity (oral) | : Not classified (Based on available data, the classification criteria are not met) |
| Acute toxicity (dermal) | : Not classified (Based on available data, the classification criteria are not met) |
| Acute toxicity (inhalation) | : Not classified (Based on available data, the classification criteria are not met) |
| Skin corrosion/irritation | : Not classified (Based on available data, the classification criteria are not met) pH: 7.2 - 7.5 (Ph. Eur. 2.2.3) |
| Serious eye damage/irritation | : Not classified (Based on available data, the classification criteria are not met) pH: 7.2 - 7.5 (Ph. Eur. 2.2.3) |
| Respiratory or skin sensitisation | : Not classified (Based on available data, the classification criteria are not met) |
| Germ cell mutagenicity | : Not classified (Based on available data, the classification criteria are not met) |
| Carcinogenicity | : Not classified (Based on available data, the classification criteria are not met) |
| Reproductive toxicity | : Not classified (Based on available data, the classification criteria are not met) |
| STOT-single exposure | : Not classified (Based on available data, the classification criteria are not met) |
| STOT-repeated exposure | : Not classified (Based on available data, the classification criteria are not met) |
| Aspiration hazard | : Not classified (Based on available data, the classification criteria are not met) |

SECTION 12: Ecological information

12.1. Toxicity

| | |
|--------------------------|---|
| Ecology - general | : Ecological problems are not known or expected under normal use. |
| Acute aquatic toxicity | : Not classified |
| Chronic aquatic toxicity | : Not classified |

12.2. Persistence and degradability

| GMP Dendritic Cell Medium (DC), Serum-free, w/o Phenol Red | |
|---|---------------------------|
| Persistence and degradability | Bioaccumulation unlikely. |

12.3. Bioaccumulative potential

| GMP Dendritic Cell Medium (DC), Serum-free, w/o Phenol Red | |
|---|----------------|
| Log Pow | Not determined |
| Log Kow | Not determined |

12.4. Mobility in soil

No additional information available

12.5. Results of PBT and vPvB assessment

| GMP Dendritic Cell Medium (DC), Serum-free, w/o Phenol Red | |
|--|--|
| This substance/mixture does not meet the PBT criteria of REACH regulation, annex XIII | |
| This substance/mixture does not meet the vPvB criteria of REACH regulation, annex XIII | |

12.6. Other adverse effects

No additional information available

SECTION 13: Disposal considerations

13.1. Waste treatment methods

| | |
|---------------------------------|---|
| Regional legislation (waste) | : This material and its container must be disposed of in a safe way, and as per local legislation. |
| Sewage disposal recommendations | : Do not allow to enter drains or water courses. |
| Additional information | : Empty the packaging completely prior to disposal. Handle uncleaned empty containers as full ones. |

SECTION 14: Transport information

In accordance with ADR / RID / IMDG / IATA / ADN

| ADR | IMDG | IATA | ADN | RID |
|---|----------------|----------------|----------------|----------------|
| 14.1. UN number | | | | |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| 14.2. UN proper shipping name | | | | |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| 14.3. Transport hazard class(es) | | | | |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| 14.4. Packing group | | | | |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| 14.5. Environmental hazards | | | | |
| Not applicable | Not applicable | Not applicable | Not applicable | Not applicable |
| No supplementary information available | | | | |

14.6. Special precautions for user

- Overland transport

Not applicable

- Transport by sea

Not applicable

- Air transport

Not applicable

- Inland waterway transport

Not applicable

- Rail transport

Not applicable

14.7. Transport in bulk according to Annex II of Marpol and the IBC Code

IBC code : Not applicable.

SECTION 15: Regulatory information

15.1. Safety, health and environmental regulations/legislation specific for the substance or mixture

15.1.1. EU-Regulations

Contains no REACH substances with Annex XVII restrictions

Contains no substance on the REACH candidate list

Contains no REACH Annex XIV substances

Other information, restriction and prohibition regulations : Export and import of dangerous chemicals: Not subject to Regulation (EC) No 649/2012. Ozone layer depleting substances: Not subject to Regulation (EC) No 1005/2009. Persistent organic pollutants (POPs): Not subject to Regulation (EC) No 850/2004. Control of major-accident hazards (COMAH, Seveso III): Not subject to Directive 2012/18/EC.

15.1.2. National regulations

No additional information available

15.2. Chemical safety assessment

Not specifically applicable
Mixtures

SECTION 16: Other information

Other information

: For this mixture, a safety data sheet is legally not required, neither due to REACH art. 31(1) nor due to REACH art. 31(3). Therefore, the compilation of this safety data sheet is done voluntarily and the safety data sheet doesn't have to fit the requirements of REACH annex II to the full extent.

This information is based on our current knowledge and is intended to describe the product for the purposes of health, safety and environmental requirements only. It should not therefore be construed as guaranteeing any specific property of the product.