4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

SECTION 1: Identification of the substance/mixture and of the company/undertaking

1.1 Product identifier

Trade name	: 4Cell® Nutri-T
Product code	: 05-11F2001-1K
Substance name	: 4Cell® Nutri-T Medium - 05-11F2001-1

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

: 4Cell® Nutri-T medium is a serum- and xeno-free medium, optimized for human T cell activation and expansion towards ex-vivo cell culture applications.
ex-vivo cell culture applications.

1.3 Details of the supplier of the safety data sheet

Company	:	Biological Industries Israel Beit-Haemek Itd. Kibbutz Beit-Haemek 2511500, Israel
Telephone	:	+972.4.9960595
Telefax	:	+972.4.9968896
E-mail address of person responsible for the SDS	:	Department CA-FM Environment@sartorius.com
Contact person	:	Product Management
Technical Support:	:	South America, Asia/Pacific/Europe/Middle East/Africa: support@bioind.com

SAFETY DATA SHEET according to Regulation (EC) No. 1907/2006

SVIFCTFAS

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

1.4 Emergency telephone number

UK/Europe/Global: +44.1865.407.333 (Sartorius 29003) United States: +1.866.928.0789 (Toll free), (Sartorius 29003) Canada: +1.800.579.7421 (Toll free), (Sartorius 29003) United States & Canada (geographic): +1.202.464.2554 (Sartorius 29003) Asia-Pacific/East & South East Asia (without China): +65.3158.1074 (Sartorius 29003) Japan: +81.3.4578.9341 (Sartorius 29003) Indonesia: 007.803.011.0293 (Toll free Indonesia), (Sartorius 29003) Malaysia: +60.3.6207.4347 (Sartorius 29003) Thailand: 001.800.120.666.751 (Toll free Thailand), (Sartorius 29003) India: 000.800.100.7479 (Toll free India) or +65.3158.1198 (Sartorius 29003) South Korea: +82.2.3479.8401 (Sartorius 29003) Australia: +61.2.8014.4558 (Sartorius 29003) New Zealand: +64.9.929.1483 (Sartorius 29003) Middle East/Africa: +44.1235.239671 (Sartorius 29003)

SECTION 2: Hazards identification

2.1 Classification of the substance or mixture

Classification (REGULATION (EC) No 1272/2008)

Not a hazardous substance or mixture according to Regulation (EC) No. 1272/2008.

2.2 Label elements

Labelling REGULATION (EC) No 1272/2008

Not a hazardous substance or mixture according to Regulation (EC) No. 1272/2008.

2.3 Other hazards

None known.

SECTION 3: Composition/information on ingredients

3.2 Mixtures

Chemical nature

: Aqueous solution Reagent

Hazardous components

Remarks

: No dangerous ingredients according to Regulation (EC) No. 1907/2006

SVIFCTSVS

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

SECTION 4: First aid measures

4.1 Description of first aid measures

If inhaled :		ve to fresh air. If irritation or breathing problems persist, ase contact medical personnel.
In case of skin contact :		sh off with plenty of water. ymptoms persist, call a physician.
In case of eye contact :		se thoroughly with plenty of water, also under the eyelids. ye irritation persists, consult a specialist.
If swallowed :		proughly rinse mouth with water. After rinsing, drink large ounts of water.
	lf s	ymptoms persist, call a physician.

4.2 Most important symptoms and effects, both acute and delayed

None known.

4.3 Indication of any immediate medical attention and special treatment needed

SECTION 5: Firefighting measures

5.1 Extinguishing media Suitable extinguishing media	: Use water spray, alcohol-resistant foam, dry chemical or car- bon dioxide.
5.2 Special hazards arising from t	the substance or mixture
Specific hazards during fire- fighting	: No information available.
5.3 Advice for firefighters	
Special protective equipment for firefighters	: In the event of fire, wear self-contained breathing apparatus.
Further information	: The product itself does not burn. Use extinguishing measures that are appropriate to local cir- cumstances and the surrounding environment.

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

SECTION 6: Accidental release measures

6.1 Personal precautions, protective equipment and emergency procedures		
Personal precautions :	Avoid formation of aerosol.	
6.2 Environmental precautions		
Environmental precautions :	Do not let product enter drains. Should not be released into the environment.	
6.3 Methods and material for contain	nment and cleaning up	
Methods for cleaning up :	Soak up with inert absorbent material (e.g. sand, silica gel, acid binder, universal binder, sawdust). Pick up and transfer to properly labelled containers.	

6.4 Reference to other sections

SECTION 7: Handling and storage

7.1 Precautions for safe handling

Advice on safe handling	: Avoid formation of aerosol. Clean contaminated surfaces thoroughly.
Hygiene measures	: Avoid contact with skin, eyes and clothing. Keep away from food, drink and animal feeds. When using do not smoke.

7.2 Conditions for safe storage, including any incompatibilities

Requirements for storage areas and containers	: Keep container tightly closed in a dry, cool (2-8 °C), dark and well-ventilated place. Protect from exposure to light.
Storage class (TRGS 510)	: 10-13, German Storage Class 10 to 13
7.3 Specific end use(s) Specific use(s)	: No data available

SECTION 8: Exposure controls/personal protection

8.1 Control parameters

8.2 Exposure controls

Personal protective equipment

Eye protection : Goggles

4Cell® Nutri-T

Versio 1.2	on	Revision Date: 25.08.2020		DS Number: AG-R0002224	Date of last issue: 20.08.2020 Date of first issue: 20.08.2020
Hand protection Remarks		:	breakthrough tim gloves. Also take tions under which	ne instructions regarding permeability and e which are provided by the supplier of the into consideration the specific local condi- n the product is used, such as the danger of nd the contact time.	
S	Skin ar	nd body protection	:	Lab coat	

SECTION 9: Physical and chemical properties

9.1 Information on basic physical and chemical properties

Appearance	: Aqueous solution
Colour	: red
Odour	: not significant
Melting point/freezing point	: 0 °C
Boiling point/boiling range	: ≥ 100 °C
Vapour pressure	: 23 hPa (20 °C)
Density	: 1 g/cm3 (20 °C)
Donoky	. i g/ointo (20 0)
Solubility(ies)	
Water solubility	: completely miscible
Viscosity Viscosity, dynamic	: 0,952 mPa.s (20 °C)
	, , , , , , , , , , , , , , , , , , ,

9.2 Other information

No data available

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

SECTION 10: Stability and reactivity

10.1 Reactivity

10.2 Chemical stability

10.3 Possibility of hazardous reactions

Hazardous reactions : None known.

10.4 Conditions to avoid

10.5 Incompatible materials

Materials to avoid

: Oxidizing agents Acids and bases

10.6 Hazardous decomposition products

No decomposition if stored and applied as directed.

SECTION 11: Toxicological information

11.1 Information on toxicological effects

Further information

Product:

Remarks: The product contains no substances known to be hazardous to health in concentrations which need to be taken into account.

SECTION 12: Ecological information

12.1 Toxicity

No data available

- **12.2 Persistence and degradability** No data available
- 12.3 Bioaccumulative potential

No data available

12.4 Mobility in soil

No data available

12.5 Results of PBT and vPvB assessment Not relevant

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

12.6 Other adverse effects

Product:

Additional ecological infor-	:	Remarks: In large quantities, this product may affect the func-
mation		tioning of waste-water treatment plants.

SECTION 13: Disposal considerations

13.1 Waste treatment methods

Product

: In accordance with local and national regulations.

SECTION 14: Transport information

14.1 UN number

14.2 UN proper shipping name

- 14.3 Transport hazard class(es)
- 14.4 Packing group
- 14.5 Environmental hazards

14.6 Special precautions for user

Remarks

: Not dangerous goods in the meaning of ADR/RID, ADN, IMDG-Code, ICAO/IATA-DGR

14.7 Transport in bulk according to Annex II of MARPOL 73/78 and the IBC Code

Not applicable for product as supplied.

SECTION 15: Regulatory information

15.1 Safety, health and environmental regulations/legislation specific for the substance or mix- ture				
Water contaminating class (Germany)	: WGK 1 slightly hazardous to water			
Volatile organic compounds	: Volatile organic compounds (VOC) content: 0,01 %			

15.2 Chemical safety assessment

SECTION 16: Other information

SAFETY DATA SHEET according to Regulation (EC) No. 1907/2006

SVIFCTFA3

4Cell® Nutri-T

Version	Revision Date:	SDS Number:	Date of last issue: 20.08.2020
1.2	25.08.2020	SAG-R0002224	Date of first issue: 20.08.2020

The information provided in this Safety Data Sheet is correct to the best of our knowledge, information and belief at the date of its publication. The information given is designed only as a guidance for safe handling, use, processing, storage, transportation, disposal and release and is not to be considered a warranty or quality specification. The information relates only to the specific material designated and may not be valid for such material used in combination with any other materials or in any process, unless specified in the text.

REG_WORLD / EN