

Page 1/6

Safety data sheet

according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

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

- · Product identifier
- · Trade name NutriCoat® Solution
- · Article number: 05-760-1-15
- *Relevant identified uses of the substance or mixture and uses advised against No further relevant information available.*
- \cdot Application of the substance / the mixture

The NutriCoat Attachment Solution is a xeno-free, ready-to-use, composed of sterile, affinity purified human fibrinogen (hFN) that supports the attachment and spreading of animal cells cultured in serum-free medium. NutriCoat Attachment Solution application validated and quality tested to provide optimal attachment and growth of animal cells, including hMSC from various sources such as bone marrow (BM-MSC), adipose tissue (AT-MSC), and umbilical cord tissue (UC-MSC).

 \cdot Details of the supplier of the safety data sheet

• Manufacturer/Supplier: Biological Industries Israel Beit Haemek Ltd. BEIT HAEMEK 25115 ISRAEL michalb@bioind.com

- · Further information obtainable from: Product safety department.
- Emergency telephone number: During normal opening times: +972/4/9962467

2 Hazards identification

· Classification of the substance or mixture

- Classification according to Regulation (EC) No 1272/2008 The product is not classified, according to the CLP regulation.
- · Label elements
- · Labelling according to Regulation (EC) No 1272/2008 Void
- · Hazard pictograms Void
- · Signal word Void
- · Hazard statements Void
- · Other hazards
- · Results of PBT and vPvB assessment
- **PBT:** Not applicable.
- · vPvB: Not applicable.

3 Composition/information on ingredients

- · Chemical characterisation: Mixtures
- · Description: Mixture of substances listed below with nonhazardous additions.
- · Dangerous components: Void
- Additional information:

All other ingredients in the proportions used are considered non-hazardous by regulating bodies worldwide

4 First aid measures

- · Description of first aid measures
- · General information: No special measures required.
- · After inhalation: Supply fresh air; consult doctor in case of complaints.

(Contd. on page 2)

GB



Page 2/6

(Contd. of page 1)

Safety data sheet according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

Trade name: NutriCoat® Solution

- After skin contact: Generally the product does not irritate the skin.
- After eye contact: Rinse opened eye for several minutes under running water.
- After swallowing: Do not induce vomiting; call for medical help immediately.
- Information for doctor:
- · Most important symptoms and effects, both acute and delayed No further relevant information available.
- · Indication of any immediate medical attention and special treatment needed
- No further relevant information available.

5 Firefighting measures

- · Extinguishing media
- Suitable extinguishing agents: Use fire extinguishing methods suitable to surrounding conditions.
- Special hazards arising from the substance or mixture No further relevant information available.
- · Advice for firefighters
- · Protective equipment: No special measures required.

6 Accidental release measures

- · Personal precautions, protective equipment and emergency procedures Not required
- · Environmental precautions: No special measures required.
- Methods and material for containment and cleaning up: No special measures required.
- · Reference to other sections
- See Section 7 for information on safe handling.
- See Section 8 for information on personal protection equipment.
- See Section 13 for disposal information.

7 Handling and storage

- · Handling:
- · Precautions for safe handling No special measures required.
- Information about fire and explosion protection: No special measures required.
- · Conditions for safe storage, including any incompatibilities
- · Storage:
- \cdot Requirements to be met by storerooms and receptacles: 15-30 °C
- · Information about storage in one common storage facility: Not Required.
- Further information about storage conditions: Not required.
- *Specific end use(s) No further relevant information available.*

8 Exposure controls/personal protection

• Additional information about design of technical facilities: No further data; see item 7.

- · Control parameters
- · Ingredients with limit values that require monitoring at the workplace:

The product does not contain any relevant quantities of materials with critical values that have to be monitored at the workplace.

• Additional information: The lists valid during the making were used as basis.



Page 3/6

Safety data sheet according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

Trade name: NutriCoat® Solution

| (Contd. of page 2) |
|---|
| Exposure controls |
| Personal protective equipment: |
| General protective and hygienic measures: |
| The usual precautionary measures are to be adhered to when handling chemicals. |
| Protection of hands: |
| The glove material has to be impermeable and ressistante to the product/substace/preperation |
| Due to missing tests no recommendation to the glove material can be given for the product/ the preparation/ |
| the chemical mixture. |
| Selection of the glove material on consideration of the penetration times, rates of diffusion and the |
| degradation |
| |

· Material of gloves

The selection of the suitable gloves does not only depend on the material, but also on further marks of quality and varies from manufacturer to manufacturer. As the product is a preparation of several substances, the resistance of the glove material can not be calculated in advance and has therefore to be checked prior to the application.

· Penetration time of glove material

The exact break through time has to be found out by the manufacturer of the protective gloves and has to be observed.

· Eye protection: Not required.

| Information on basic physical and chem | ical properties | |
|--|---|--|
| General Information | | |
| Appearance: | | |
| Form: | Liquid | |
| Colour: | Colorless | |
| Odour: | Characteristic | |
| Odour threshold: | Not determined. | |
| pH-value: | Not applicable. | |
| Change in condition | | |
| Melting point/freezing point: | Undetermined. | |
| Initial boiling point and boiling range: | Undetermined. | |
| Flash point: | Not applicable. | |
| Flammability (solid, gas): | Not applicable. | |
| Decomposition temperature: | Not determined. | |
| Auto-ignition temperature: | Product is not selfigniting. | |
| Explosive properties: | Product does not present an explosion hazard. | |
| Explosion limits: | | |
| Lower: | Not determined. | |
| Upper: | Not determined. | |
| Vapour pressure: | Not determined. | |
| Density: | Not determined. | |
| Relative density | Not determined. | |
| Vapour density | Not determined. | |

GB



Page 4/6

Safety data sheet according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

Trade name: NutriCoat® Solution

| | | (Contd. of page 3 |
|---|--|-------------------|
| · Evaporation rate | Not determined. | |
| · Solubility in / Miscibility with | | |
| water: | Fully miscible. | |
| · Partition coefficient: n-octanol/water: | Not determined. | |
| · Viscosity: | | |
| Dynamic: | Not determined. | |
| Kinematic: | Not determined. | |
| · Solvent content: | | |
| Water: | >95.0 % | |
| VOC (EC) | 0.00 % | |
| • Other information | No further relevant information available. | |

10 Stability and reactivity

· Reactivity No further relevant information available.

- · Chemical stability
- Thermal decomposition / conditions to be avoided: No decomposition if used according to specifications.
- Possibility of hazardous reactions No dangerous reactions known.
- · Conditions to avoid No further relevant information available.
- *Incompatible materials:* No further relevant information available.
- · Hazardous decomposition products: No dangerous decomposition products known.

11 Toxicological information

· Information on toxicological effects

- · Acute toxicity Based on available data, the classification criteria are not met.
- · Primary irritant effect:
- · Skin corrosion/irritation Based on available data, the classification criteria are not met.
- Serious eye damage/irritation Based on available data, the classification criteria are not met.
- Respiratory or skin sensitisation Based on available data, the classification criteria are not met.
- · CMR effects (carcinogenity, mutagenicity and toxicity for reproduction)
- Germ cell mutagenicity Based on available data, the classification criteria are not met.
- · Carcinogenicity Based on available data, the classification criteria are not met.
- Reproductive toxicity Based on available data, the classification criteria are not met.
- STOT-single exposure Based on available data, the classification criteria are not met.
- STOT-repeated exposure Based on available data, the classification criteria are not met.
- · Aspiration hazard Based on available data, the classification criteria are not met.

12 Ecological information

· Toxicity

- Aquatic toxicity: No further relevant information available.
- · Persistence and degradability No further relevant information available.
- · Behaviour in environmental systems:
- · Bioaccumulative potential No further relevant information available.
- · Mobility in soil No further relevant information available.

(Contd. on page 5)

GB -



Page 5/6

(Contd. of page 4)

Safety data sheet according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

Trade name: NutriCoat® Solution

· Additional ecological information:

· General notes: Generally not hazardous for water

· Results of PBT and vPvB assessment

- **PBT:** Not applicable.
- · vPvB: Not applicable.

· Other adverse effects No further relevant information available.

13 Disposal considerations

· Waste treatment methods

· Recommendation Smaller quantities can be disposed of with household waste.

· Uncleaned packaging:

· Recommendation: Disposal must be made according to official regulations.

| UN-Number | not reculated | |
|---|-----------------|--|
| ADR, ADN, IMDG, IATA | not regulated | |
| UN proper shipping name ADR, ADN, IMDG, IATA | not regulated | |
| Transport hazard class(es) | | |
| ADR, ADN, IMDG, IATA | | |
| Class | not regulated | |
| Packing group | | |
| ADR, IMDG, IATA | not regulated | |
| Environmental hazards: | Not applicable. | |
| Special precautions for user | Not applicable. | |
| Transport in bulk according to Annex I | I of Marpol | |
| and the IBC Code | Not applicable. | |

15 Regulatory information

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

· Directive 2012/18/EU

· Named dangerous substances - ANNEX I None of the ingredients is listed.

· Chemical safety assessment: A Chemical Safety Assessment has not been carried out.

16 Other information

This information is based on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship.

· Department issuing SDS: Product safety department.

(Contd. on page 6)

GB



Page 6/6

Safety data sheet according to 1907/2006/EC, Article 31

Printing date 14.11.2018

Version 1

Revision date: 12.11.2018

Trade name: NutriCoat® Solution

(Contd. of page 5)

• Contact: Ms. BADICHI GITLIN • Abbreviations and acronyms:

ADR: Accord européen sur le transport des marchandises dangereuses par Route (European Agreement concerning the International Carriage of Dangerous Goods by Road) IMDG: International Maritime Code for Dangerous Goods

IATA: International Air Transport Association

GHS: Globally Harmonised System of Classification and Labelling of Chemicals

EINECS: European Inventory of Existing Commercial Chemical Substances

ELINCS: European List of Notified Chemical Substances

CAS: Chemical Abstracts Service (division of the American Chemical Society)

VOC: Volatile Organic Compounds (USA, EU)

PBT: Persistent, Bioaccumulative and Toxic vPvB: very Persistent and very Bioaccumulative