

## **Safety data sheet**

**according to 1907/2006/EC, Article 31**

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

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

· **Product identifier**

· **Trade name:** MSC NutriStem® XF Supplement Mix

· **Article number:** 05-201-1

· **Relevant identified uses of the substance or mixture and uses advised against**  
No further relevant information available.

· **Application of the substance / the mixture**

MSC NutriStem® XF Medium is a defined, serum-free, xeno-free culture medium designed for optimal growth and expansion of human mesenchymal stem/stromal cells (hMSC) derived from a variety of sources, including bone marrow (BM-MSC), adipose tissue (AT-MSC) and umbilical cord matrix (UC-MSC). MSC NutriStem® XF Medium supports long-term growth of hMSC while maintaining their self-renewal and multi-lineage differentiation potential.

· **Details of the supplier of the safety data sheet**

· **Manufacturer/Supplier:**

Biological Industries Israel Beit Haemek Ltd.  
BEIT HAEMEK 2511500  
ISRAEL  
Michal.badihi@sartorius.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)

## **Safety data sheet**

**according to 1907/2006/EC, Article 31**

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

**Trade name: MSC NutriStem® XF Supplement Mix**

(Contd. of page 1)

- **After skin contact:** Generally the product does not irritate the skin.
- **After eye contact:**  
Rinse opened eye for several minutes under running water. If symptoms persist, consult a doctor.
- **After swallowing:** Wash mouth with water if person is conscious and seek medical advice.
- **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 protective equipment required.

### **6 Accidental release measures**

- **Personal precautions, protective equipment and emergency procedures**  
Use personal protective equipment. Clean contaminated surfaces thoroughly.
- **Environmental precautions:** Use personal protective equipment. Clean contaminated surfaces thoroughly.
- **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 precautions are necessary if used correctly.  
Avoid freeze/thaw cycles
- **Information about fire - and explosion protection:** No special measures required.
- **Conditions for safe storage, including any incompatibilities**
- **Storage:**
- **Requirements to be met by storerooms and receptacles:** (-10)-(-20) °C
- **Information about storage in one common storage facility:** Not required.
- **Further information about storage conditions:** Avoid exposure to light.
- **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.

(Contd. on page 3)

GB

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

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

**Trade name: MSC NutriStem® XF Supplement Mix**

(Contd. of page 2)

- **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.*
- **Exposure controls**
- **Personal protective equipment:**
- **General protective and hygienic measures:**  
*The usual precautionary measures are to be adhered to when handling chemicals.*
- **Respiratory protection:** *Not required*
- **Protection of hands:**  
*The glove material has to be impermeable and resistant to the product/substance/preparation  
 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:** *Safety glasses*

## 9 Physical and chemical properties

- **Information on basic physical and chemical properties**
- **General Information**
- **Appearance:**

<b>Form:</b>	Solution
<b>Colour:</b>	Clear
<b>Odour:</b>	Uncharacteristic.
<b>Odour threshold:</b>	Not determined.
- **pH-value:** *Not applicable.*
- **Change in condition**

<b>Melting point/freezing point:</b>	Undetermined.
<b>Initial boiling point and boiling range:</b>	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:**

<b>Lower:</b>	Not determined.
---------------	-----------------

(Contd. on page 4)

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

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

**Trade name: MSC NutriStem® XF Supplement Mix**

(Contd. of page 3)

<b>Upper:</b>	<i>Not determined.</i>
· <b>Vapour pressure at 20 °C:</b>	23 hPa
· <b>Density:</b>	<i>Not determined.</i>
· <b>Relative density</b>	<i>Not determined.</i>
· <b>Vapour density</b>	<i>Not determined.</i>
· <b>Evaporation rate</b>	<i>Not determined.</i>
· <b>Solubility in / Miscibility with water:</b>	Fully miscible.
· <b>Partition coefficient: n-octanol/water:</b>	<i>Not determined.</i>
· <b>Viscosity:</b>	
<b>Dynamic:</b>	<i>Not determined.</i>
<b>Kinematic:</b>	<i>Not determined.</i>
· <b>Solvent content:</b>	
<b>Water:</b>	>70.0 %
<b>VOC (EC)</b>	0.00 %
· <b>Solids content:</b>	<1.0 %
· <b>Other information</b>	<i>No further relevant information available.</i>

### 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 (carcinogenicity, 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.*

GB

(Contd. on page 5)

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

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

**Trade name: MSC NutriStem® XF Supplement Mix**

(Contd. of page 4)

### 12 Ecological information

- **Toxicity**
- **Aquatic toxicity:** Not toxic to the aquatic life
- **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.
- **Additional ecological information:**
- **General notes:**  
 Water hazard class 1 (German Regulation) (Self-assessment): slightly hazardous for water  
 Do not allow undiluted product or large quantities of it to reach ground water, water course or sewage system.
- **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** No special recommendation
- **Uncleaned packaging:**
- **Recommendation:** Disposal must be made according to official regulations.
- **Recommended cleansing agents:** Water, if necessary together with cleansing agents.

### 14 Transport information

- |   |                 |
|---|-----------------|
| · <b>UN-Number</b>  |                 |
| · <b>ADR, ADN, IMDG, IATA</b>   | not regulated   |
| · <b>UN proper shipping name</b>  |                 |
| · <b>ADR, ADN, IMDG, IATA</b>   | not regulated   |
| · <b>Transport hazard class(es)</b>   |                 |
| · <b>ADR, ADN, IMDG, IATA</b>   |                 |
| · <b>Class</b>  | not regulated   |
| · <b>Packing group</b>  |                 |
| · <b>ADR, IMDG, IATA</b>  | not regulated   |
| · <b>Environmental hazards:</b>   |                 |
| · <b>Marine pollutant:</b>  | No              |
| · <b>Special precautions for user</b>                                       | Not applicable. |
| · <b>Transport in bulk according to Annex II of Marpol and the IBC Code</b> | Not applicable. |

(Contd. on page 6)

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

Printing date 12.07.2020

Version 2

Revision date: 22.11.2017

**Trade name: MSC NutriStem® XF Supplement Mix**

(Contd. of page 5)

· **UN "Model Regulation":** not regulated

**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.
- **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*