

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

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

1.1 Product identifier

Trade name : Mix PBS
Substance name : Mix PBS

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

1.3 Details of the supplier of the safety data sheet

Company : Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Göttingen

E-mail address of person responsible for the SDS : Department CA-FM Environment@sartorius.com

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 dangerous substance according to GHS.

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

2.3 Other hazards

None known.

SECTION 3: Composition/information on ingredients

3.2 Mixtures

Chemical nature : Laboratory chemicals
Solid
Mixture of sodium chloride

Hazardous components

SECTION 4: First aid measures

4.1 Description of first aid measures

If inhaled : Fresh Air. If disease continues, please contact a medical expert.

In case of skin contact : If on skin, rinse well with water.

In case of eye contact : Rinse immediately with plenty of water, also under the eyelids. If eye irritation persists, consult a specialist.

If swallowed : Do NOT induce vomiting.
Call a physician immediately.
Rinse mouth with water.
Give small amounts of water to drink.

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 : Water
Foam
Carbon dioxide (CO₂)
Dry powder

5.2 Special hazards arising from the substance or mixture

Specific hazards during fire-fighting : Hazardous gas or vapour can develop in case of fire.

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

5.3 Advice for firefighters

Special protective equipment for firefighters : Wear self-contained breathing apparatus for firefighting if necessary.

Further information : Standard procedure for chemical fires.
Use extinguishing measures that are appropriate to local circumstances and the surrounding environment.

SECTION 6: Accidental release measures

6.1 Personal precautions, protective equipment and emergency procedures

6.2 Environmental precautions

Environmental precautions : Do not let product enter drains.

6.3 Methods and material for containment and cleaning up

Methods for cleaning up : Sweep up and shovel into suitable container for disposal.

6.4 Reference to other sections

SECTION 7: Handling and storage

7.1 Precautions for safe handling

Hygiene measures : Do not smoke. When using do not eat or drink.

7.2 Conditions for safe storage, including any incompatibilities

Requirements for storage areas and containers : Keep container tightly closed in a dry and well-ventilated place.

7.3 Specific end use(s)

SECTION 8: Exposure controls/personal protection

8.1 Control parameters

8.2 Exposure controls

Personal protective equipment

Eye protection : Safety glasses

Hand protection

Remarks : Nitrile rubber

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

Skin and body protection : Lab coat

SECTION 9: Physical and chemical properties

9.1 Information on basic physical and chemical properties

Appearance : solid, crystalline
Colour : white
Odour : odourless
pH : ca. 6
Vapour pressure : 1.013 hPa
Density : 2,17 g/cm³ (20 °C)

9.2 Other information

No data available

SECTION 10: Stability and reactivity

10.1 Reactivity

10.2 Chemical stability

10.3 Possibility of hazardous reactions

Hazardous reactions : No dangerous reaction known under conditions of normal use.

10.4 Conditions to avoid

Conditions to avoid : Heat, flames and sparks.

10.5 Incompatible materials

Materials to avoid : Oxidizing agents

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

10.6 Hazardous decomposition products

SECTION 11: Toxicological information

11.1 Information on toxicological effects

Further information

Product:

Remarks: The product contains no substances which at their given concentration, are considered to be hazardous to health.

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

12.6 Other adverse effects

Product:

Additional ecological information : Remarks: This product has no known ecotoxicological effects.

SECTION 13: Disposal considerations

13.1 Waste treatment methods

Product : In accordance with local and national regulations.

Mix PBS

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	27.11.2019	SAG-R0002069	Date of first issue: 27.11.2019

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 mixture

Water contaminating class (Germany) : WGK 1 slightly hazardous to water

15.2 Chemical safety assessment

SECTION 16: Other information

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.

DE / EN