



SAFETY DATA SHEET

1.0 IDENTIFICATION

Product Name	Streptavidin-HRP
Manufacturer	Quanterix Corporation
Address	900 Middlesex Turnpike, Building 1, Billerica, MA 01821
Product Number	1858263
Telephone	(617) 301-9400
Intended Use of the Product	
Immunoassay on Quanterix SP-X; For Research Use Only. Not for Diagnostic Purposes.	

2.0 HAZARD(S) IDENTIFICATION

2.1 Classification

While this material is not considered hazardous by the OSHA Hazard Communication Standard (29 CFR 1910.1200), this SDS contains valuable information critical to the safe handling and proper use of the product. This SDS should be retained and available for employees and other users of this product.

2.2 GHS Label Elements

Not applicable.

2.3 Signal Word

Not applicable.

2.4 Hazard Statements

No known significant effects or critical hazards.

2.5 Precautionary Statements

May be harmful by inhalation, ingestion, or skin adsorption. May cause eye, skin, or respiratory system irritation. To the best of our knowledge, the toxicological properties have not been thoroughly investigated.

2.6 Hazards Not Otherwise Classified

Not applicable.



3.0 COMPOSITION – INFORMATION ON INGREDIENTS

3.1 Substance / Mixture

Mixture.

3.2 Ingredient Name^{1,2}

1. Streptavidin-HRP

COMPONENT	PERCENTAGE	CAS #
Peroxidase-conjugated Streptavidin	-	N/A
Sodium Phosphate	-	7558-79-4
Sodium Chloride	-	7647-14-5
Bovine Serum Albumin	-	9048-46-8
Non-Hazardous Proprietary Ingredient	-	N/A

NOTE 1: N/A means not applicable or proprietary information. The specific chemical identity and/or exact percentage of composition have been withheld as a trade secret [29CFR1910.1200].

NOTE 2: There are no additional ingredients present which, within the current knowledge of the supplier and in the concentrations applicable, are classified as hazardous to health or the environment and hence require reporting in this section.

4.0 FIRST AID MEASURES

4.1 Description of Necessary First Aid Measures

1. General

Never give anything by mouth to an unconscious person. If you feel unwell, seek medical advice (show the label where possible).

2. Eye Contact

Immediately flush eyes with plenty of water, occasionally lifting the upper and lower eyelids. Check for and remove any contact lenses. Get medical attention if irritation occurs.

3. Inhalation

Remove victim to fresh air and keep at rest in a position comfortable for breathing. Get medical attention if symptoms occur.

4. Skin Contact

Flush contaminated skin with plenty of water. Remove contaminated clothing and shoes. Get medical attention if symptoms occur.

5. Ingestion

Wash out mouth with water. Remove victim to fresh air and keep at rest in a position comfortable for breathing. If material has been swallowed and the exposed person is conscious, give small



quantities of water to drink. Do not induce vomiting unless directed to do so by medical personnel. Get medical attention if symptoms occur

4.2 Most Important Symptoms and Effects, Acute and Delayed

1. Potential Acute Health Effects

Eye Contact: No known significant effects or critical hazards.

Inhalation: No known significant effects or critical hazards.

Skin Contact: No known significant effects or critical hazards.

Ingestion: No known significant effects or critical hazards.

2. Over-exposure Signs and Symptoms

Eye Contact: No data available.

Inhalation: No data available.

Skin Contact: No data available.

Ingestion: No data available.

5.0 FIRE-FIGHTING MEASURES

5.1 Extinguishing Media

1. Suitable Extinguishing Media

Use an extinguishing agent suitable for the surrounding fire.

2. Unsuitable Extinguishing Media

None known.

5.2 Special Hazards Arising from the Substance or Mixture

1. Fire Hazard

Not flammable.

2. Explosion Hazard

Product is not explosive.

3. Thermal Decomposition Products

Decomposition products may include the following materials:

- carbon dioxide;
- carbon monoxide;
- sulfur oxides;
- metal oxide/oxides



5.3 Advice for Fire-Fighters

1. Special Protective Actions

Promptly isolate the scene by removing all persons from the vicinity of the incident if there is a fire. No action shall be taken involving any personal risk or without suitable training.

2. Special Protective Equipment

Fire-fighters should wear appropriate protective equipment and self-contained breathing apparatus (SCBA) with a full face-piece operated in positive pressure mode.

6.0 ACCIDENTAL RELEASE MEASURES

6.1 Personal Precautions, Protective Equipment and Emergency Procedures

1. For Non-Emergency Personnel

No action shall be taken involving any personal risk or without suitable training. Evacuate surrounding areas. Keep unnecessary and unprotected personnel from entering. Do not touch or walk through spilled material. Put on appropriate personal protective equipment.

2. For Emergency Responders

If specialized clothing is required to deal with the spillage, take note of any information in Section 8 on suitable and unsuitable materials. See also the information in "For nonemergency personnel".

6.2 Environmental Precautions

Avoid dispersal of spilled material and runoff and contact with soil, waterways, drains and sewers. Inform the relevant authorities if the product has caused environmental pollution (sewers, waterways, soil or air).

6.3 Methods and Materials for Containment and Cleaning Up

Stop leak if without risk. Move containers from spill area. Dilute with water and mop up if water-soluble. Alternatively, or if water-insoluble, absorb with an inert dry material and place in an appropriate waste disposal container. Dispose of via a licensed waste disposal contractor.

7.0 HANDLING AND STORAGE

7.1 Precautions for Safe Handling

Put on appropriate personal protective equipment (see section 8). Eating, drinking and smoking should be prohibited in areas where this material is handled, stored and processed. Workers should wash hands and face before eating, drinking and smoking. Do not ingest. Avoid contact with eyes, skin and clothing. Keep in the original container or an approved alternative made from a compatible material, kept tightly closed when not in use. Empty containers retain product residue and can be hazardous. Do not reuse container.



7.2 Conditions for Safe Storage, Including Incompatibilities

Store in accordance with local regulations. Store in original container protected from direct sunlight in a dry, cool and well-ventilated area, away from incompatible materials (see section 10) and food and drink. Keep container tightly closed and sealed until ready for use. Containers that have been opened must be carefully resealed and kept upright to prevent leakage. Do not store in unlabeled containers. Use appropriate containment to avoid environmental contamination.

8.0 **EXPOSURE CONTROLS AND PERSONAL PROTECTION**

8.1 Occupational Exposure Limits

Ingredient Name	Exposure Limits
n/a	n/a

8.2 Recommended Monitoring procedures:

If this product contains ingredients with exposure limits, personal, workplace atmosphere or biological monitoring may be required to determine the effectiveness of the ventilation or other control measures and/or the necessity to use respiratory protective equipment.

8.3 Engineering Measures:

No special ventilation requirements. Good general ventilation should be sufficient to control worker exposure to airborne contaminants. If this product contains ingredients with exposure limits, use process enclosures, local exhaust ventilation or other engineering controls to keep worker exposure below any recommended or statutory limits.

8.4 Hygiene Measures

Wash hands, forearms and face thoroughly after handling chemical products, before eating, smoking and using the lavatory and at the end of the working period. Appropriate techniques should be used to remove potentially contaminated clothing. Wash contaminated clothing before reusing. Ensure that eyewash stations and safety showers are close to the workstation location.

8.5 Personal Protection

1. **Respiratory**

Use a properly fitted, air-purifying or air-fed respirator complying with an approved standard if a risk assessment indicates this is necessary. Respirator selection must be based on known or anticipated exposure levels, the hazards of the product and the safe working limits of the selected respirator.

2. **Hands**

Chemical-resistant, impervious gloves complying with an approved standard should be worn at all times when handling chemical products if a risk assessment indicates this is necessary.



3. Eyes

Safety eyewear complying with an approved standard should be used when a risk assessment indicates this is necessary to avoid exposure to liquid splashes, mists, gases or dusts.

4. Skin

Personal protective equipment for the body should be selected based on the task being performed and the risks involved and should be approved by a specialist before handling this product.

9.0 PHYSICAL AND CHEMICAL PROPERTIES

9.1 Information on Basic Physical and Chemical Properties

1. Appearance

Physical State	Solid
Color	Light Brown
Odor	Odorless
Odor Threshold	No data available.
pH	7.6 when rehydrated with indicated volume of H ₂ O
Melting Point	No data available.
Boiling Point	No data available.
Flash Point	No data available.
Evaporation Rate	No data available.
Flammability (solid, gas)	No data available.
Lower and Upper Explosive Limits	No data available.
Vapor Pressure	No data available.
Vapor Density	No data available.
Relative Density	No data available.
Solubility	No data available.
Partition Coefficient: n-octanol/water	No data available.
Auto-ignition Temperature	No data available.
Decomposition Temperature	No data available.



Viscosity

No data available.

10.0 STABILITY AND REACTIVITY

10.1 Reactivity

No specific test data related to reactivity available for this product or its ingredients.

10.2 Chemical Stability

The product is stable.

10.3 Possibility of Hazardous Reactions

Under normal conditions of storage and use, hazardous reactions will not occur.

10.4 Conditions to Avoid

Keep away from heat.

10.5 Incompatible Materials

None known.

10.6 Hazardous Decomposition Products

Under normal conditions of storage and use, hazardous decomposition products should not be produced.

11.0 TOXICOLOGICAL INFORMATION

11.1 Information on Toxicological Effects

1. Acute Toxicity

No data available.

2. Irritation / Corrosion

No data available.

3. Sensitization

No data available.

4. Mutagenicity

No data available.

5. Carcinogenicity

No data available.

6. Reproductive Toxicity

No data available.



7. Teratogenicity

No data available.

8. Specific Organ Toxicity (Single Exposure)

No data available.

9. Specific Organ Toxicity (Repeated Exposure)

No data available.

10. Aspiration Hazard

No data available.

11. Information on the Likely Routes of Exposure

No data available.

12. Potential Acute Health Effects

<u>Eye Contact:</u>	No known significant effects or critical hazards.
<u>Inhalation:</u>	No known significant effects or critical hazards.
<u>Skin Contact:</u>	No known significant effects or critical hazards.
<u>Ingestion:</u>	No known significant effects or critical hazards.

13. Symptoms Related to the Physical, Chemical and Toxicological Characteristics

<u>Eye Contact:</u>	No known significant effects or critical hazards.
<u>Inhalation:</u>	No known significant effects or critical hazards.
<u>Skin Contact:</u>	No known significant effects or critical hazards.
<u>Ingestion:</u>	No known significant effects or critical hazards.

14. Potential Chronic Health Effects

<u>General</u>	No known significant effects or critical hazards.
<u>Carcinogenicity:</u>	No known significant effects or critical hazards.
<u>Mutagenicity</u>	No known significant effects or critical hazards.
<u>Teratogenicity:</u>	No known significant effects or critical hazards.
<u>Developmental:</u>	No known significant effects or critical hazards.
<u>Fertility Effects</u>	No known significant effects or critical hazards.

12.0 ECOLOGICAL INFORMATION

12.1 Ecological Toxicity

Avoid release into environment. Runoff from fire control or dilution water may cause pollution.



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 Other Adverse Effects

No data available.

13.0 DISPOSAL CONSIDERATIONS

13.1 Waste Treatment Methods

The generation of waste should be avoided or minimized wherever possible. Dispose of surplus and non-recyclable products via a licensed waste disposal contractor. Disposal of this product, solutions and any by-products should at all times comply with the requirements of environmental protection and waste disposal legislation and any regional local authority requirements. Avoid dispersal of spilled material and runoff and contact with soil, waterways, drains and sewers.

14.0 TRANSPORTATION INFORMATION

14.1 Hazard Class/ UN Number/ Proper Shipping Name

	ADR/RID	ADN/ADNR	IMDG	IATA
UN Number	Not regulated for transport	Not regulated for transport	Not regulated for transport	Not regulated for transport
UN proper shipping name	Not regulated for transport	Not regulated for transport	-	-
Transport hazard class(es)	Not regulated for transport	Not regulated for transport	-	-
Packing group	-	-	-	-
Environmental hazards	No	No	No	No
Special precaution for user	Not regulated for transport	Not regulated for transport	Not regulated for transport	Not regulated for transport

14.2 DOT

Transport in bulk according to Annex II of MARPOL 73/78 and the IBC Code: Not available

15.0 REGULATORY INFORMATION

No data available.

16.0 OTHER INFORMATION

The information and recommendations contained herein are based upon tests believed to be reliable. However, Quanterix Corporation does not guarantee their accuracy or completeness NOR SHALL ANY OF THIS INFORMATION CONSTITUTE A WARRANTY, WHETHER EXPRESSED OR

IMPLIED, AS TO THE SAFETY OF THE GOODS, THE MERCHANTABILITY OF THE GOODS, OR THE FITNESS OF THE GOODS FOR A PARTICULAR PURPOSE. Adjustment to conform to actual conditions of usage may be required. Quanterix Corporation assumes no responsibility for results obtained or for incidental or consequential damages, including lost profits arising from the use of these data. No warranty against infringement of any patent, copyright or trademark is made or implied.