

# **SAFETY DATA SHEET**

## **1.0 IDENTIFICATION**

<b>Product Name</b>	<b>SARS-CoV-2 Spike IgG Advantage Kit</b>
<b>Manufacturer</b>	Quanterix Corporation
<b>Address</b>	900 Middlesex Turnpike, Building 1, Billerica, MA 01821
<b>Product Number</b>	<b>103769</b>
<b>Telephone</b>	(617) 301-9400
<b>Intended Use of the Product</b> For Research Use Only. Not for Diagnostic Purposes.	

## **2.0 HAZARDS IDENTIFICATION**

### **Classification of the Substance or Mixture**

#### **Classification (GHS-US)**

Not classified

### **Label Elements**

#### **GHS-US Labeling**

Not applicable

### **Other Hazards**

No additional information available

### **Unknown Acute Toxicity (GHS-US)**

No data available



## 3.0 COMPOSITION – INFORMATION ON INGREDIENTS<sup>1</sup>

### Bead Reagent

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	90-100%	Not classified
Tris(hydroxymethyl)aminomethane	77-86-1	0.1-1%	201-064-4
Tris(hydroxymethyl)aminomethane hydrochloride	118-53-1	0.1-1%	214-684-5
Sodium Chloride	7647-14-5	0.1-1%	231-598-3
Ethylenediaminetetraacetic acid disodium salt dihydrate	6381-92-6	0.1-1%	200-449-4
Bovine Serum Albumin	9048-46-8	1-10%	232-936-2
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.1-1%	Unlisted
Tween 20	9005-64-5	0.1-1%	Unlisted
Non-Hazardous Proprietary Ingredient	N/A	-	Unlisted

### Detector Reagent

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	90-100%	Not classified
Sodium phosphate, dibasic	7558-79-4	1-10%	231-448-7
Potassium phosphate, Monobasic	7778-77-0	0.1-1%	231-913-4
Sodium Chloride	7647-14-5	0.5-5.0%	231-598-3
Potassium chloride	7447-40-7	0.01-0.1%	231-211-8
Ethylenediaminetetraacetic acid disodium salt dihydrate	6381-92-6	0.1-1%	200-449-4
Bovine Serum Albumin	9048-46-8	1-10%	232-936-2
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.01-0.1%	Unlisted
Tween 20	9005-64-5	0.1-1.0%	Unlisted
Non-Hazardous Proprietary Ingredient	N/A	–	Unlisted

## Sample Diluent

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	66-90%	Not classified
Sodium phosphate, dibasic	7558-79-4	1-5%	231-448-7
Potassium phosphate, Monobasic	7778-77-0	1-5%	231-913-4
Potassium chloride	7447-40-7	1-5%	231-211-8
Sodium Chloride	7647-14-5	1-5%	231-598-3
Fetal Bovine Serum	Unlisted	1-10%	Unlisted
Triton X-100	9002-93-1	0.01-0.1%	Unlisted
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.01-0.1%	Unlisted
Bovine Serum Albumin	9048-46-8	1-10%	232-936-2
Nonylphenol, ethoxylated	9016-45-9	0.1-1%	500-024-6

## SBG Reagent

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	90-100%	Not classified
Sodium phosphate, dibasic	7558-79-4	1-5%	231-448-7
Potassium phosphate, Monobasic	7778-77-0	0.5-2%	231-913-4
Potassium chloride	7447-40-7	0.1 – 1%	231-211-8
Sodium Chloride	7647-14-5	0.1-1%	231-598-3
Ethylenediaminetetraacetic acid disodium salt dihydrate	6381-92-6	0.1-1%	200-449-4
Bovine Serum Albumin	9048-46-8	1-10%	232-936-2
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.1-0.5%	Unlisted
Tween 20	9005-64-5	0.1-1%	Unlisted
Magnesium chloride	7786-30-3	0.1-1%	Unlisted
Non-Hazardous Proprietary Ingredient	N/A	-	Unlisted

## RGP Reagent

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	90-100%	Not classified
Sodium phosphate, dibasic	7558-79-4	0.1-0.5%	231-448-7
Potassium phosphate, Monobasic	7778-77-0	0.01-0.1%	231-913-4
Potassium chloride	7447-40-7	0.01-0.1%	231-211-8
Sodium Chloride	7647-14-5	0.1-1%	231-598-3
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.1-0.5%	Unlisted
Pluronic F-127	9003-11-6	0.01-0.1%	Unlisted
Resorufin b-galactopyranoside	95079-19-9	0.001-0.01%	Unlisted

## Calibrators & Controls

Component	CAS No.	Percent Composition	EINECS number
Water	7732-18-5	66-90%	Not classified
Sodium phosphate, dibasic	7558-79-4	1-5%	231-448-7
Potassium phosphate, Monobasic	7778-77-0	1-5%	231-913-4
Potassium chloride	7447-40-7	1-5%	231-211-8
Sodium Chloride	7647-14-5	1-5%	231-598-3
Fetal Bovine Serum	Unlisted	1-10%	Unlisted
Triton X-100	9002-93-1	0.01-0.1%	Unlisted
Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone	55965-84-9	0.01-0.1%	Unlisted
Bovine Serum Albumin	9048-46-8	1-10%	232-936-2
Nonylphenol, ethoxylated	9016-45-9	0.1-1%	500-024-6
Non-Hazardous Proprietary Ingredient	N/A	-	Unlisted

**NOTE 1:** These ingredients, within the current knowledge of the supplier and in the concentrations applicable, are not classified as hazardous to health or to the environment.



## 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

<b>Physical State</b>	No data available.
<b>Color</b>	No data available.
<b>Odor</b>	No data available.
<b>Odor Threshold</b>	No data available.
<b>pH</b>	No data available.
<b>Melting Point</b>	No data available.
<b>Boiling Point</b>	No data available.
<b>Flash Point</b>	No data available.
<b>Evaporation Rate</b>	No data available.
<b>Flammability (solid, gas)</b>	No data available.
<b>Lower and Upper Explosive Limits</b>	No data available.
<b>Vapor Pressure</b>	No data available.
<b>Vapor Density</b>	No data available.
<b>Relative Density</b>	No data available.
<b>Solubility</b>	No data available.
<b>Partition Coefficient: n- octanol/water</b>	No data available.
<b>Auto-ignition Temperature</b>	No data available.
<b>Decomposition Temperature</b>	No data available.
<b>Viscosity</b>	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**

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 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**

Not regulated.

### **14.2 DOT**

Not regulated.

### **14.3 IATA**

Not regulated.

### **14.4 IMDG**

Not regulated.

## **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.