

# Safety Data Sheet

## 1. CHEMICAL PRODUCT AND COMPANY IDENTIFICATION

|   |   |
|---|---|
| <b>Product Name</b>                                 | <b>Homebrew Detector Diluent</b>  |
| <b>Description</b>                                  | A water-based buffer to be used on the Quanterix instrument (Simoa) for the detection and quantification of purified proteins. Kit consists of a mixture of reagents that are non-hazardous, non-toxic and non-infectious compounds; not for human use. |
| <b>Manufacturer</b>                                 | Quanterix Corporation   |
| <b>Address</b>                                      | 900 Middlesex Turnpike, Building 1, Billerica, MA 01821   |
| <b>Product Number</b>                               | <b>101363</b>   |
| <b>SDS Issued</b>                                   | 20 May 2019   |
| <b>Telephone</b>                                    | (617) 301-9400  |
| <b>Intended Use of the Product</b>                  |   |
| For Research Use Only. Not for Diagnostic Purposes. |   |

## 2. 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. COMPOSITION – INFORMATION ON INGREDIENTS**

**Homebrew Detector/Sample Diluent**

| Component  | CAS No.    | Percent Composition | EINECS number  |
|--|------------|---------------------|----------------|
| Water  | 7732-18-5  | 90-100%             | Not classified |
| Sodium phosphate, dibasic  | 7558-79-4  | 0.2-2%              | 231-448-7      |
| Potassium phosphate, Monobasic   | 7778-77-0  | 0.03-0.3%           | 231-913-4      |
| Sodium Chloride  | 7647-14-5  | 0.5-5%              | 231-598-3      |
| Potassium chloride   | 7447-40-7  | 0.01-0.1%           | 231-211-8      |
| Bovine Serum Albumin   | 9048-46-8  | 0.05-0.5%           | 232-936-2      |
| Tween 20   | 9005-64-5  | 0.1-1%              | Unlisted       |
| Ethylenediaminetetraacetic acid disodium salt dihydrate                            | 6381-92-6  | 0.1-1%              | 200-449-4      |
| Interference blockers  | N/A        | N/A                 | Unlisted       |
| Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone | 55965-84-9 | 0.05-0.5%           | Unlisted       |

**4. FIRST AID MEASURES**

**Description of First Aid Measures**

**First-aid Measures General:** Never give anything by mouth to an unconscious person. If you feel unwell, seek medical advice(show the label where possible).

**First-aid Measures After Inhalation:** When symptoms occur: go into open air and ventilate suspected area.

**First-aid Measures After Skin Contact:** Remove contaminated clothing. Drench affected area with water for at least 15 minutes.

**First-aid Measures After Eye Contact:** Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

**First-aid Measures After Ingestion:** Rinse mouth. Do NOT induce vomiting.

**Most important symptoms and effects, both acute and delayed**

**Symptoms/Injuries:** None expected under normal conditions of use.

**Symptoms/Injuries After Skin Contact:** Contact during a long period may cause light irritation.

**Symptoms/Injuries After Eye Contact:** Direct contact with the eyes is likely irritating.

**Symptoms/Injuries After Ingestion:** May be harmful if swallowed.

**Indication of Any Immediate Medical Attention and Special Treatment Needed**

If medical advice is needed, have product container or label at hand.

## 5. FIRE FIGHTING MEASURES

---

### 5.1. Extinguishing Media

**Suitable Extinguishing Media:** Alcohol foam, polymer foam, dry chemical powder, carbon dioxide, water spray, fog.

**Unsuitable Extinguishing Media:** None known.

### 5.2. Special Hazards Arising From the Substance or Mixture

**Fire Hazard:** Not flammable.

**Explosion Hazard:** Product is not explosive.

**Reactivity:** If the product is involved in a fire, it can release toxic chlorine gases.

### 5.3. Advice for Firefighters

**Firefighting Instructions:** Exercise caution when fighting any chemical fire.

**Protection During Firefighting:** Do not enter fire area without proper protective equipment, including respiratory protection.

## 6. ACCIDENTAL RELEASE MEASURES

---

### 6.1. Personal Precautions, Protective Equipment and Emergency Procedures

**General Measures:** Do not breathe fumes from fires or vapours from decomposition.

#### 6.1.1. For Non-emergency Personnel

**Protective Equipment:** Use appropriate personal protection equipment (PPE).

**Emergency Procedures:** Evacuate unnecessary personnel.

#### 6.1.2. For Emergency Responders

**Protective Equipment:** Equip cleanup crew with proper protection.

**Emergency Procedures:** Ventilate area.

### 6.2. Environmental Precautions

Prevent entry to sewers and public waters.

### 6.3. Methods and Material for Containment and Cleaning Up

**For Containment:** Absorb and/or contain spill with inert material, then place in suitable container.

**Methods for Cleaning Up:** Clear up spills immediately and dispose of waste safely.

### 6.4. Reference to Other Sections

See heading 8, Exposure Controls and Personal Protection.

## 7. HANDLING AND STORAGE

**Handling:** Refer to section 8. Wear appropriate personal protective equipment when using this product.

**Storage:** Storage at 2-8 degrees Celsius

## 8. EXPOSURE CONTROLS, PERSONAL PROTECTION

### 8.1. Control Parameters

No Occupational Exposure Limits (OELs) have been established for this product or its chemical components.

### 8.2. Exposure Controls

#### Appropriate Engineering Controls

: Emergency eye wash fountains should be available in the immediate vicinity of any potential exposure.

#### Personal Protective Equipment

: Gloves. In case of splash hazard: safety glasses.



#### Hand Protection

: Wear chemically resistant protective gloves.

#### Eye Protection

: Chemical goggles or safety glasses.

#### Other Information

: When using, do not eat, drink or smoke.

## 9. PHYSICAL AND CHEMICAL PROPERTIES

|                   |                   |
|-------------------|-------------------|
| Appearance:       | Clear Liquid      |
| Odor:             | Odorless          |
| Solubility:       | Miscible in water |
| Physical state:   | Liquid            |
| Melting Point:    | Not Available     |
| Boiling Point:    | Not Available     |
| Vapor Pressure:   | Not Available     |
| Vapor Density:    | Not Available     |
| pH:               | 7.0-8.0           |
| Evaporation Rate: | Not Available     |

## 10. STABILITY AND REACTIVITY

|                                       |            |
|---------------------------------------|------------|
| Incompatibility: (materials to avoid) | None known |
| Stability:                            | Stable     |
| Hazardous Polymerization:             | None       |
| Conditions to Avoid:                  | None       |

## 11. TOXICOLOGICAL INFORMATION

### 11.1. Information On Toxicological Effects

Acute Toxicity : Not classified

|  |  |
|--|--|
| <b>Potassium chloride (7447-40-7)</b>  |  |
| LD50 Oral Rat  | 2600 mg/kg                                 |
| <b>Dihydrogen potassium phosphate (7778-77-0)</b>  |  |
| LD50 Dermal Rabbit   | > 4640 mg/kg                               |
| <b>Sodium chloride (7647-14-5)</b>   |  |
| LD50 Oral Rat  | 3 g/kg                                     |
| LD50 Dermal Rabbit   | > 10 g/kg                                  |
| LC50 Inhalation Rat (mg/l)   | > 42 g/m <sup>3</sup> (Exposure time: 1 h) |
| <b>Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone (55965-84-9)</b> |  |
| LD50 Oral Rat  | 53 mg/kg                                   |

Skin Corrosion/Irritation: Not classified

Serious Eye Damage/Irritation: Not classified

Respiratory or Skin Sensitization: Not classified

Germ Cell Mutagenicity: Not classified

Carcinogenicity: Not classified

Reproductive Toxicity: Not classified

Specific Target Organ Toxicity (Single Exposure): Not classified

Specific Target Organ Toxicity (Repeated Exposure): Not classified

Aspiration Hazard: Not classified

Symptoms/Injuries After Skin Contact: Contact during a long period may cause light irritation.

Symptoms/Injuries After Eye Contact: Direct contact with the eyes is likely irritating.

Symptoms/Injuries After Ingestion: May be harmful if swallowed.

## 12. ECOLOGICAL INFORMATION

Data not yet available.

## 13. DISPOSAL CONSIDERATIONS

### 13.1. Waste treatment methods

**Waste Disposal Recommendations:** Dispose of waste material in accordance with all local, regional, national and international regulations.

## 14. TRANSPORTATION INFORMATION

To the best of our knowledge, this product is not regulated as a hazardous material.

## 15. REGULATORY INFORMATION

#### US Federal Regulations

|   |
|---|
| <b>Potassium chloride (7447-40-7)</b>                                     |
| Listed on the United States TSCA (Toxic Substances Control Act) inventory |
| <b>Dihydrogen potassium phosphate (7778-77-0)</b>                         |
| Listed on the United States TSCA (Toxic Substances Control Act) inventory |
| <b>Sodium chloride (7647-14-5)</b>  |
| Listed on the United States TSCA (Toxic Substances Control Act) inventory |
| <b>Water (7732-18-5)</b>  |
| Listed on the United States TSCA (Toxic Substances Control Act) inventory |
| <b>Polyoxyethylene sorbitan monolaurate (9005-64-5)</b>                   |
| Listed on the United States TSCA (Toxic Substances Control Act) inventory |

#### US State Regulations

|   |
|---|
| <b>Potassium chloride (7447-40-7)</b>                   |
| U.S. - Texas - Effects Screening Levels - Long Term     |
| U.S. - Texas - Effects Screening Levels - Short Term    |
| <b>Sodium chloride (7647-14-5)</b>                      |
| U.S. - Texas - Effects Screening Levels - Long Term     |
| U.S. - Texas - Effects Screening Levels - Short Term    |
| <b>Polyoxyethylene sorbitan monolaurate (9005-64-5)</b> |
| U.S. - Texas - Effects Screening Levels - Long Term     |
| U.S. - Texas - Effects Screening Levels - Short Term    |

## 16. OTHER INFORMATION

### Disclaimer:

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.