

# Simoa Wash Buffer B

Quanterix

## Safety Data Sheet

According to Federal Register / Vol. 77, No. 58 / Monday, March 26, 2012 / Rules and Regulations; Revision Date: 20th May, 2019

Version: 1.0

### SECTION 1: IDENTIFICATION OF THE SUBSTANCE/MIXTURE AND OF THE COMPANY

#### 1.1. Product Identifier

**Product Form:** Mixtures

**Product Name:** Simoa Wash Buffer B

**Item/Product Number:** Item #102867 (Single unit) and 103079 (2 Pack)

#### 1.2. Intended Use of the Product

**Use of the Substance/Mixture:** For use with Quanterix system. For Research Use Only. Not for Diagnostic Purposes.

#### 1.3. Name, Address, and Telephone of the Responsible Party

##### Company

Quanterix Corporation  
900 Middlesex Turnpike, Building 1,  
Billerica, MA 01821  
T 617-301-9400  
[www.quanterix.com](http://www.quanterix.com)

### SECTION 2: HAZARDS IDENTIFICATION

#### 2.1. Classification of the Substance or Mixture

##### Classification (GHS-US)

Not classified

#### 2.2. Label Elements

##### GHS-US Labeling

Not applicable

#### 2.3. Other Hazards

No additional information available

#### 2.4. Unknown Acute Toxicity (GHS-US):

No data available

### SECTION 3: COMPOSITION/INFORMATION ON INGREDIENTS

#### 3.1. Mixtures

| Name   | Product Identifier  | %        | Classification (GHS-US)   |
|--|---------------------|----------|---|
| Water  | (CAS No) 7732-18-5  | 90 - 100 | Not classified  |
| Sodium chloride  | (CAS No) 7647-14-5  | 1 - 5    | Not classified  |
| Disodium orthophosphate heptahydrate   | (CAS No) 7782-85-6  | 1 - 5    | Skin Irrit. 2, H315   |
| Dihydrogen potassium phosphate   | (CAS No) 7778-77-0  | 0.01-1   | Eye Irrit. 2A, H319   |
| Saccharide*  | (CAS No) 57-50-1    | 3-30     | Not classified  |
| Potassium chloride   | (CAS No) 7447-40-7  | 0.01 - 1 | Eye Irrit. 2B, H320   |
| Mixture, 3(2H)-isothiazolone, 5-chloro-2-methyl- with 2-methyl-3(2H)-isothiazolone | (CAS No) 55965-84-9 | 0.01 - 1 | Acute Tox. 3 (Oral), H301<br>Acute Tox. 3 (Dermal), H311<br>Acute Tox. 3 (Inhalation), H331<br>Skin Corr. 1B, H314<br>Skin Sens. 1, H317<br>Aquatic Chronic 1, H410 |
| Surfactant*  | (CAS No) 9005-64-5  | 0.01 - 1 | Not classified  |

\*SPECIFICS OF THE COMPOUND ARE QUANTERIX TRADEMARK INFORMATION AND CANNOT BE DISCLOSED.

Full text of H-phrases: see section 16

## SECTION 4: FIRST AID MEASURES

### 4.1. 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.

### 4.2. 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.

### 4.3. Indication of Any Immediate Medical Attention and Special Treatment Needed

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

## SECTION 5: FIREFIGHTING 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.

## SECTION 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.

## SECTION 7: HANDLING AND STORAGE

### 7.1. Precautions for Safe Handling

**Hygiene Measures:** Handle in accordance with good industrial hygiene and safety procedures. Wash hands and other exposed areas with mild soap and water before eating, drinking, or smoking and again when leaving work.

### 7.2. Conditions for Safe Storage, Including Any Incompatibilities

**Storage Conditions:** Store in a dry, cool and well-ventilated place. Keep container closed when not in use.

**Incompatible Products:** Strong acids. Strong bases. Strong oxidizers.

## 7.3. Specific End Use(s)

Immunoassay on Simoa HD-1.

## SECTION 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.

## SECTION 9: PHYSICAL AND CHEMICAL PROPERTIES

### 9.1. Information on Basic Physical and Chemical Properties

|  |                     |
|--|---------------------|
| Physical State                             | : Liquid            |
| Appearance                                 | : Clear. Colorless. |
| Odor                                       | : No data available |
| Odor Threshold                             | : No data available |
| pH   | : No data available |
| Relative Evaporation Rate (butylacetate=1) | : No data available |
| Melting Point                              | : No data available |
| Freezing Point                             | : No data available |
| Boiling Point                              | : No data available |
| Flash Point                                | : No data available |
| Auto-ignition Temperature                  | : No data available |
| Decomposition Temperature                  | : No data available |
| Flammability (solid, gas)                  | : No data available |
| Vapor Pressure                             | : No data available |
| Relative Vapor Density at 20 °C            | : No data available |
| Relative Density                           | : No data available |
| Specific Gravity                           | : Not available     |
| Solubility                                 | : No data available |
| Log Pow                                    | : No data available |
| Log Kow                                    | : No data available |
| Viscosity, Kinematic                       | : No data available |
| Viscosity, Dynamic                         | : No data available |
| Explosive Properties                       | : No data available |
| Oxidizing Properties                       | : No data available |
| Explosive Limits                           | : Not applicable    |

### 9.2. Other Information

No additional information available

## SECTION 10: STABILITY AND REACTIVITY

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

**Chemical Stability:** Product is stable.

**Possibility of Hazardous Reactions:** Hazardous polymerization will not occur.

**Conditions to Avoid:** Direct sunlight. Extremely high or low temperatures.

**Incompatible Materials:** Strong acids. Strong bases. Strong oxidizers.

**Hazardous Decomposition Products:** Carbon oxides (CO, CO<sub>2</sub>). Upon heating, toxic fumes are formed. (chlorine)

**SECTION 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.

**SECTION 12: ECOLOGICAL INFORMATION****12.1. Toxicity**

|                                       |   |
|---------------------------------------|---|
| <b>Potassium chloride (7447-40-7)</b> |   |
| LC50 Fish 1                           | 1060 mg/l (Exposure time: 96 h - Species: Lepomis macrochirus [static])                     |
| EC50 Daphnia 1                        | 825 mg/l (Exposure time: 48 h - Species: Daphnia magna)                                     |
| EC50 Other Aquatic Organisms 1        | 2500 mg/l (Exposure time: 72 h - Species: Desmodemus subspicatus)                           |
| LC 50 Fish 2                          | 750 - 1020 mg/l (Exposure time: 96 h - Species: Pimephales promelas [static])               |
| EC50 Daphnia 2                        | 83 mg/l (Exposure time: 48 h - Species: Daphnia magna [Static])                             |
| <b>Sodium chloride (7647-14-5)</b>    |   |
| LC50 Fish 1                           | 5560 (5560 - 6080) mg/l (Exposure time: 96 h - Species: Lepomis macrochirus [flow-through]) |
| EC50 Daphnia 1                        | 1000 mg/l (Exposure time: 48 h - Species: Daphnia magna)                                    |
| LC 50 Fish 2                          | 12946 mg/l (Exposure time: 96 h - Species: Lepomis macrochirus [static])                    |
| EC50 Daphnia 2                        | 340.7 (340.7 - 469.2) mg/l (Exposure time: 48 h - Species: Daphnia magna [Static])          |

**12.2. Persistence and Degradability**

|                               |                  |
|-------------------------------|------------------|
| <b>System Wash Buffer 2</b>   |                  |
| Persistence and Degradability | Not established. |

**12.3. Bioaccumulative Potential**

|                                    |                      |
|------------------------------------|----------------------|
| <b>System Wash Buffer 2</b>        |                      |
| Bioaccumulative Potential          | Not established.     |
| <b>Sodium chloride (7647-14-5)</b> |                      |
| BCF fish 1                         | (no bioaccumulation) |

**12.4. Mobility in Soil** No additional information available

# Wash Buffer B

## Safety Data Sheet

according to Federal Register / Vol. 77, No. 58 / Monday, March 26, 2012 / Rules and Regulations

### 12.5. Other Adverse Effects

**Other Information** : Avoid release to the environment.

## SECTION 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.

## SECTION 14: TRANSPORT INFORMATION

In Accordance With ICAO/IATA/DOT/TDG

**14.1. UN Number** Not applicable

**14.2. UN Proper Shipping Name** Not regulated for transport.

### 14.3. Additional Information

**Other information** : No supplementary information available.

**Transport by Sea** Not regulated for transport.

**Air Transport** Not regulated for transport.

## SECTION 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    |

## SECTION 16: OTHER INFORMATION

**Other Information** : This document has been prepared in accordance with the SDS requirements of the OSHA Hazard Communication Standard 29 CFR 1910.1200.

### GHS Full Text Phrases:

|                           |  |
|---------------------------|--|
| Acute Tox. 3 (Dermal)     | Acute toxicity (dermal) Category 3                               |
| Acute Tox. 3 (Inhalation) | Acute toxicity (inhalation) Category 3                           |
| Acute Tox. 3 (Oral)       | Acute toxicity (oral) Category 3                                 |
| Aquatic Chronic 1         | Hazardous to the aquatic environment - Chronic Hazard Category 1 |
| Eye Irrit. 2A             | Serious eye damage/eye irritation Category 2A                    |
| Eye Irrit. 2B             | Serious eye damage/eye irritation Category 2B                    |
| Skin Corr. 1B             | Skin corrosion/irritation Category 1B                            |
| Skin Irrit. 2             | Skin corrosion/irritation Category 2                             |

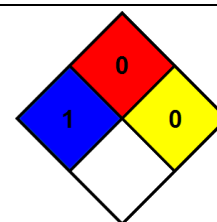
# Wash Buffer B

## Safety Data Sheet

according to Federal Register / Vol. 77, No. 58 / Monday, March 26, 2012 / Rules and Regulations

|              |   |
|--------------|---|
| Skin Sens. 1 | Skin sensitization Category 1                               |
| STOT SE 3    | Specific target organ toxicity (single exposure) Category 3 |
| H301         | Toxic if swallowed  |
| H311         | Toxic in contact with skin                                  |
| H314         | Causes severe skin burns and eye damage                     |
| H315         | Causes skin irritation                                      |
| H317         | May cause an allergic skin reaction                         |
| H319         | Causes serious eye irritation                               |
| H320         | Causes eye irritation                                       |
| H331         | Toxic if inhaled  |
| H335         | May cause respiratory irritation                            |
| H410         | Very toxic to aquatic life with long lasting effects        |

- NFPA Health Hazard** : 1 - Exposure could cause irritation but only minor residual injury even if no treatment is given.
- NFPA Fire Hazard** : 0 - Materials that will not burn.
- NFPA Reactivity** : 0 - Normally stable, even under fire exposure conditions, and are not reactive with water.



### HMIS III Rating

- Health** : 1 Slight Hazard - Irritation or minor reversible injury possible
- Flammability** : 0 Minimal Hazard
- Physical** : 0 Minimal Hazard

*This information is based on our current knowledge and is intended to describe the product for the purposes of health, safety and environmental requirements only. It should not therefore be construed as guaranteeing any specific property of the product*

SDS US (GHS HazCom) - US Only