

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

**1.1. Product Identifier**

**Product Name:** SR-X Sealing Oil; Item number 102767

**1.2. Intended Use of the Product**

**Use of the substance/mixture:** Consumable used for Immunoassay testing; For Research Use Only. Not for Diagnostic Purposes.

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

**Company**

Quanterix Corporation  
113 Hartwell Avenue  
Lexington, MA 02421  
T 617-301-9400  
[www.quanterix.com](http://www.quanterix.com)

**1.4. Emergency Telephone Number**

For Chemical Emergency, Spill, Leak, Fire, Exposure, or Accident, call CHEMTREC – Day or Night (800) 424-9300

**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):**

Not applicable

**SECTION 3: COMPOSITION/INFORMATION ON INGREDIENTS**

**3.1. Substances**

Name	Product Identifier	%	Classification (GHS-US)
Fluorinated Oil*	Cannot be disclosed*	100	Not classified

\* Specifics of this compound is Quanterix Trademark and cannot be disclosed, but all the relevant safety information is 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 Inhalation:** Not expected to present a significant inhalation hazard under anticipated conditions of normal use.

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

**Symptoms/Injuries After Eye Contact:** May cause slight irritation to eyes.

# System Sealing Oil

## Safety Data Sheet

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

**Symptoms/Injuries After Ingestion:** May cause gastrointestinal irritation, nausea, vomiting and diarrhoea.

### 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. Use extinguishing media appropriate for surrounding fire.

**Unsuitable Extinguishing Media:** A heavy water stream may spread burning liquid. Do not use water jet to extinguish.

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

**Fire Hazard:** Not flammable.

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

**Reactivity:** Depolymerization may occur in the presence of some metal oxides at temperatures above 288°C (550°F).

### 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:** No special measures required.

#### 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 testing

## 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**

: Use ventilation when the oil is heated above 260°C (550°F).

**Personal Protective Equipment**

: Gloves. In case of splash hazard: safetyglasses.



**Hand Protection**

: Wear chemically resistant protective gloves.

# System Sealing Oil

## Safety Data Sheet

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

**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** : Viscous. Clear, colorless, oily liquid.  
**Odor** : Odorless  
**Odor Threshold** : No data available  
**pH** : 7  
**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** : Does not ignite. PMCC.  
**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** : 1.86 - 1.91 @24°C  
**Specific Gravity** : Not available  
**Solubility** : Negligible.  
**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  
**Pour Point** : -45 to -5°C (-49 to -25°F)

**9.2. Other Information** No additional information available

## SECTION 10: STABILITY AND REACTIVITY

**Reactivity:** Depolymerization may occur in the presence of some metal oxides at temperatures above 288°C (550°F).

**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>). Heating above 350°C (662°F) may form potentially toxic fluorine compounds. Decomposition occurs at increasing rates as temperature is raised above 355°C (670°F).

## SECTION 11: TOXICOLOGICAL INFORMATION

### 11.1. Information On Toxicological Effects

**Acute Toxicity** : Not classified

<b>LD50 Oral Rat</b>	25000 mg/kg
<b>LD50 Dermal Rabbit</b>	17000 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

# System Sealing Oil

## Safety Data Sheet

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

**Specific Target Organ Toxicity (Repeated Exposure):** Not classified

**Aspiration Hazard:** Not classified

**Symptoms/Injuries After Inhalation:** Not expected to present a significant inhalation hazard under anticipated conditions of normal use.

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

**Symptoms/Injuries After Eye Contact:** May cause slight irritation to eyes.

**Symptoms/Injuries After Ingestion:** May cause gastrointestinal irritation, nausea, vomiting and diarrhoea.

## SECTION 12: ECOLOGICAL INFORMATION

**12.1. Toxicity:** No additional information available

**12.2. Persistence and Degradability**

System Sealing Oil	
Persistence and Degradability	Not established.

**12.3. Bioaccumulative Potential**

System Sealing Oil	
Bioaccumulative Potential	Not established.

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

**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**

Sealing Oil

Listed on the United States TSCA (Toxic Substances Control Act) inventory

**US State Regulations**

Sealing Oil

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. Not classified (Dermal)	Acute toxicity (dermal) Not classified
Acute Tox. Not classified (Oral)	Acute toxicity (oral) Not classified

# System Sealing Oil

## Safety Data Sheet

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

### NFPA Health Hazard

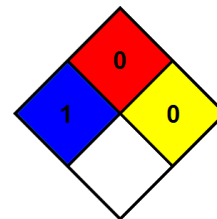
: 1 - Exposure could cause irritation but only minor residual injury even if no treatment is given.

### 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

: 0 Minimal Hazard - No significant risk to health

#### 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