

# TREVIGEN®

## MATERIAL SAFETY DATA SHEET

### PRODUCT INFORMATION

PRODUCT NAME: Phenol Red (solution)  
CATALOG NUMBER: 3430-50-01; 3430-50-06  
Chemical Name: Phenol, 4,4'- 7-Hydroxy-3H-penoxazin-3-one-10-oxide  
Physical State: Liquid in amber microvials

### EMERGENCY OVERVIEW

HMIS RATING HEALTH: 2 FLAMMABILITY: 0 PHYSICAL HAZARD: 0  
NFPA RATING HEALTH: 2 FLAMMABILITY: 0 REACTIVITY: 0

Irritant. Irritating to eyes, respiratory system and skin.

CAS#: 62758-13-8 SARA 313: No

### FIRE FIGHTING MEASURES

FLASH POINT N/A  
AUTOIGNITION TEMP N/A  
FLAMMABILITY N/A  
EXTINGUISHING MEDIA

Suitable: Water spray, Carbon dioxide, dry chemical powder, or appropriate foam.

### PROCEDURE(S) OF PERSONAL PRECAUTION(S)

Wear laboratory coat, safety goggles, rubber boots, and heavy rubber gloves.

### METHODS FOR CLEANING UP

Absorb on sand or vermiculite and place in closed containers for disposal. Ventilate area and wash spill site after material pickup is complete.

### HANDLING

Storage: Room Temperature

### STORAGE

Suitable: Keep tightly closed.

### STABILITY

Stable: Stable.  
Materials to Avoid: Strong Oxidizing agents.

### ROUTE OF EXPOSURE

Inhalation: May be harmful; material is irritating to mucous membranes and upper respiratory tract.  
Skin: May be harmful if absorbed through the skin; causes skin irritation.  
Eyes: Causes irritation.  
Ingestion: May be harmful if swallowed.

### FIRST AID

In case of contact: Immediately wash with plenty of soap and water.  
In case of eye contact: Flush eyes with water for at least 15 minutes.  
If swallowed: Never give anything by mouth to an unconscious person. Rinse mouth out with water and call a physician.

### DISPOSAL

Contact a licensed professional waste disposal service to dispose of this material. Dissolve or mix this liquid with a combustible solvent and burn in a chemical incinerator equipped with an afterburner and scrubber. Observe all federal, state, and local environmental regulations.

CONTINUED ON NEXT PAGE

Trevigen, Inc. • 8405 Helgerman Court • Gaithersburg, MD 20877 • Tel: 1-800-873-8443  
Fax: 301-560-4973 • email: info@trevigen.com

**TRANSPORT INFORMATION**

DOT: Not Hazardous  
IATA: Not Hazardous

**REGULATORY INFORMATION:****EU ADDITIONAL CLASSIFICATION:**

Symbol of Danger: Xi  
Indication of Danger: Irritant  
R: 36/37/38  
Risk Statements: Irritating to eyes, respiratory system and skin.  
S: 26  
Safety Statements: In case of contact with eyes, rinse with plenty of water and seek medical advice.

**US CLASSIFICATION AND LABEL TEXT:**

Indication of Danger: Irritant  
Risk Statements: Irritating to eyes, respiratory system and skin.  
Safety Statements: In case of contact with eyes, rinse with plenty of water and seek medical advice.  
SARA LISTED: No

**CANADA CLASSIFICATION:**

**WHMIS:** This product has been classified in accordance with the hazard criteria of the CPR, and the MSDS contains all the information required by the CPR.  
DSL: Yes  
NDSL: No

For research use only; not for drug, household, or other uses.

The above information is believed to be correct, however, the physical, chemical, and toxicological properties of the product may not yet have been fully characterized. Trevigen recommends the use of gloves, lab coats and eye protection while using or handling this product. Trevigen assumes no liability for damage resulting from handling or contact with the above product.