



<b>4</b> FIRST-AID MEASURES	<b>Includes important symptoms/ effects, acute, delayed; required treatment.</b>
<p>Skin contact: Rinse with plenty of water. If symptoms arise, call a physician.</p> <p>Eye contact: Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. If symptoms persist, call a physician.</p> <p>Ingestion: Never give anything by mouth to an unconscious person. If symptoms persist, call a physician. Do not induce vomiting without medical advice.</p> <p>Inhalation: Move to fresh air. If symptoms persist, call a physician. If not breathing, give artificial respiration.</p> <p>Notes to physician: Treat symptomatically.</p>	
<b>5</b> FIRE-FIGHTING MEASURES	<b>Lists suitable extinguishing techniques, equipment; chemical hazards from fire.</b>
<p>Suitable extinguishing media: Water spray, alcohol-resistant foam, dry chemical or CO<sub>2</sub>.</p> <p>Special protective equipment for firefighters: Wear self-contained breathing apparatus and protective suit.</p>	
<b>6</b> ACCIDENTAL RELEASE MEASURES	<b>Lists emergency procedures; protective equipment; proper methods of containment and cleanup.</b>
<p>Personal precautions: Evacuate Area. Wear self-contained breathing apparatus, rubber boots and heavy rubber gloves.</p> <p>Methods for cleaning up: Absorb with sand or vermiculite, sweep up, place in a bag and hold for waste disposal. Avoid raising dust. Ventilate area and wash spill site after material pickup is complete.</p> <p>Environmental precautions: Prevent product from entering drains.</p>	
<b>7</b> HANDLING AND STORAGE	<b>Lists precautions for safe handling and storage, including incompatibilities.</b>
<p>Handling: Should be handled by trained personnel observing good laboratory practices. Avoid breathing vapor, skin contact, and swallowing. May cause allergic reaction in sensitized individuals.</p> <p>Storage: Keep in properly labeled containers.</p>	
<b>8</b> EXPOSURE CONTROLS/ PERSONAL PROTECTION	<b>Lists OSHA's Permissible Exposure Limits (PELs); ACGIH Threshold Limit Values (TLVs); and any other exposure limit used or recommended by the chemical manufacturer, importer, or employer preparing the SDS where available as well as appropriate engineering controls; personal protective equipment (PPE).</b>
<p>Engineering measures: Ensure adequate ventilation, especially in confined areas.</p> <p>Personal protective equipment: Personal Protective Equipment requirements are dependent on the user institution's risk assessment and are specific to the risk assessment for each laboratory where this material may be used.</p> <p>Respiratory protection: In case of insufficient ventilation, wear suitable respiratory equipment.</p> <p>Hand protection: Impervious gloves.</p> <p>Eye protection: Safety glasses with side-shields.</p> <p>Skin and body protection: Lightweight protective clothing.</p> <p>Hygiene measures: Handle in accordance with good industrial hygiene and safety practice.</p>	

**9 PHYSICAL AND CHEMICAL PROPERTIES** Lists the chemical's characteristics.

**General information**

Form: No information available.

**Important health and safety environmental information**

Boiling point: No information available.

Melting point/range: No information available.

Flash point: No information available.

Autoignition temperature: No information available.

Oxidizing properties: No information available.

Water solubility: No data available.

**10 STABILITY AND REACTIVITY** Lists chemical stability and possibility of hazardous reactions.

Stability: Stable under normal conditions.

Materials to avoid: No dangerous reaction known under conditions of normal use.

Hazardous decomposition products: No information available.

Polymerization: Hazardous polymerization does not occur.

**11 TOXICOLOGICAL INFORMATION** Includes routes of exposure; related symptoms, acute and chronic effects; numerical measures of toxicity

Acute toxicity: Not hazardous.

**Principle routes of exposure/potential health effects**

Eyes: No information available.

Skin: No information available.

Inhalation: No information available.

Ingestion: May be harmful by ingestion.

Carcinogenic effects: No information available.

Mutagenic effects: No information available.

Reproductive toxicity: No information available.

Sensitization: No information available.

Target organ effects: No information available.

**12 ECOLOGICAL INFORMATION**

Ecotoxicity effects: No information available.

Mobility: No information available.

Biodegradation: Inherently biodegradable.

Bioaccumulation: Does not bioaccumulate.

**13 DISPOSAL CONSIDERATIONS** Dispose of in accordance with local regulations.

**14 TRANSPORT INFORMATION** IATA.

Proper shipping name: Not classified as dangerous in the meaning of transport regulations.  
 Hazard class: No information available.  
 Subsidiary class: No information available.  
 Packing group: No information available.  
 UN-No: No information available.

**15 REGULATORY INFORMATION** This product has been classified according to the hazard criteria of the CPR and the SDS contains all of the information required by the CPR.

**US Federal Regulations**

SARA 313: This product is not regulated by SARA.  
 Clean air act, section 112 hazardous air pollutants (HAPs) (see 40 CFR 61): This product does not contain HAPs.

**US State Regulations**

California Proposition 65: This product does not contain chemicals listed under Proposition 65.  
 WHMIS Hazard Class: Non-controlled.

**16 OTHER INFORMATION** Includes the date of preparation or last revision.

*For research use only. Not intended for human or animal diagnostic or therapeutic uses.*

The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may present unknown hazards and should be used with caution. Since the Company cannot control the actual methods, volumes, or conditions of use, the Company shall not be held liable for any damages or losses resulting

from the handling or from contact with the product as described herein. THE INFORMATION IN THIS SDS DOES NOT CONSTITUTE A WARRANTY, EXPRESSED OR IMPLIED, INCLUDING ANY IMPLIED WARRANTY OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE.

Date of preparation: November 16, 2015  
 Revision date: May 3, 2017

**END OF SAFETY DATA SHEET**