

Product: Nonidet® P-40 Substitute [Nonylphenyl-polyethylene glycol] [LU 5.11.09/MK]

<u>Components</u>	<u>Catalog No.</u>	<u>Quantity</u>
Nonidet® P-40 Substitute [Nonylphenyl-polyethylene glycol]	DG001	5 x 10ml
	DG002	10 x 10ml
	DG501	1 x 50ml
	DG514	1 x 100ml

Material Safety Data Sheet

Section 1: Chemical Identification

Name: Nonidet® P-40 Substitute [Nonylphenyl-polyethylene glycol]

Catalog #: DG001, DG002, DG501, and DG514 indicated above

Section 2: Composition/Information on Ingredients

Nonidet® P-40 Substitute [Nonylphenyl-polyethylene glycol] CAS #: 9016-45-9

Chemical Formula C₁₉H₃₂O₃

Section 3: Hazards Identification

Label precautionary statements. Irritating to respiratory system. Risk of serious damage to eyes. Possible sensitizer. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing.

Section 4: First-Aid Measures

In case of eye contact: Immediately flush eyes with copious amounts of water for at least 15 minutes.

In case of skin contact: Immediately wash skin with soap and copious amounts of water.

In case of Inhalation: Remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen.

In case of swallowing: Wash out mouth with water provided person is conscious. Call a physician.

Section 5: Fire Fighting Measures

Extinguishing Agents:

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

Special fire fighting procedures:

Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

Unusual fire and explosions hazards: Emits toxic fumes under fire conditions.

Section 6: Accidental Release Measures

Human Safety Precautions

Wear respirator, chemical safety goggles, rubber boots and heavy rubber gloves.

Cleaning/Collecting Measures:

Sweep up, place in a bag and hold for waste disposal. Ventilate area and wash spill site after material pickup is complete.

Section 7: Handling and Storage

Refer to section 8.

Section 8: Exposure Controls/Personal Protection

Personal protective equipment:

Chemical safety goggles, compatible chemical-respirator, NIOSH/MSHA-approved respirator, and safety shower and eye bath.

General Protective and Hygenic Measures:

Mechanical exhaust required. Do not breathe in dust. Avoid contact with eyes, skin and clothing. Wash thoroughly after handling. Irritant. Keep tightly closed. Store in a cool dry place.

Section 9: Physical and Chemical Properties

Physical Properties

Liquid

Section 10: Stability and Reactivity

Stability

Stable

Incompatibilities

Strong oxidizing agents

Hazardous Combustion or Decomposition Products

Carbon monoxide, carbon dioxide

Hazardous Polymerization

Not reported

Section 11: Toxicological Information

Acute Effects/Toxicity: May cause skin irritation.

Primary Irritant Effects: Harmful if absorbed through the skin, inhaled or ingested.

On the skin: Material is irritating to mucous membranes and upper respiratory tract.

On the eye: Causes eye irritation.

Toxicological properties have not been thoroughly investigated. Prolonged or repeated exposure may cause allergic reaction to certain sensitive individuals.

Additional toxicological information:

Toxicity Data

ORL - RAT LD50: 2590 UL/KG

SKN - RBT LD50: 2830 UL/KG

Only selected Registry of toxic effects of chemical substances (RTECS) data are presented here. See actual entry of RTECS for complete information.

Irritation Data

SKN-HMN 15MG/3D Remarks: Mild irritation effect

EYE-RAT 20MG Remarks: Severe irritation effects

EYE-RBT 5mg Remarks: Severe irritation effects

SKN- RBT 500mg Remarks: Mild irritation effects

EYE-GPG 20mg Remarks: Severe irritation effects

Section 12: Ecological Information

Data not yet available

Section 13: Disposal Considerations

Recommendation:

Appropriate Method of Disposal of Substance or Preparation: Must not be disposed of together with household garbage. Do not allow product to reach sewage system. Hand over to hazardous waste disposers. Contact a licensed professional waste disposal service to dispose of this material.. Dissolve or mix the material with combustible solvent and burn in a chemical incinerator equipped with an afterburner and scrubber.

Observe all federal, state and local environmental regulations.

Section 14: Transport Information

Not dangerous goods. Safe for all transport.

Section 15: Regulatory information

Irritating to eyes, respiratory system and skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing.

Section 16: Other Information

The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. *G-Biosciences/Genotech* shall not be held liable for any damage resulting from handling or from contact with the above product. See reverse side of invoice or packing slip for additional terms and conditions of sale.

G-Biosciences/Genotech St. Louis, MO, USA

Web: www.GBiosciences.com, email: technical@GBiosciences.com