

Active Motif,
1914 Palomar Oaks Way, Suite 150,
Carlsbad, CA 92008.
Phone: 760-431-1263/877-222-9543
Fax: 760-431-1351

MATERIAL SAFETY DATA SHEET (MSDS)

IDENTIFICATION:

The following products from Active Motif are covered under this MSDS.

- Purified polyclonal antibodies of mouse, rabbit or goat origin.

Components common to all of these products are buffer salts and human or animal proteins. None of these components are considered to be hazardous. The hazardous substance, sodium azide, is present in these products in very small amounts as an anti-microbial preservative.

HAZARDOUS INGREDIENTS:

<u>Materials</u>	Sodium azide, less than or equal to 0.1% (w/v) CAS# 26628-22-8
<u>Toxicity Data</u>	LD ₅₀ =27mg/kg (oral-rat)

The concentration of sodium azide is below the regulatory threshold limits according to OSHA standard 29CFR 1910.1200. However, precautions for handling potentially dangerous reagents should be practiced when using these products. To aide in determining handling procedures, we offer the following additional information.

- Physical State of sodium azide-containing antibody: Odorless, light gold color, clear liquid

The following applies to sodium azide powder

PHYSICAL DATA:

- Physical State: Odorless, white solid
- Boiling point: Undetermined
- Melting point: Undetermined
- Vapor pressure: Not applicable
- Vapor density: Not applicable
- Solubility in water: Soluble

REACTIVITY DATA:

The product has no known dangerous reactions, nor known dangerous decomposition products. Avoid strong oxidizers.

FIRE AND EXPLOSION HAZARD DATA:

There is no unusual fire or explosion hazard associated with this product.

- Flash point: Not applicable
- Flammable Limits in Air: Not applicable
- Extinguishing media: Water, carbon dioxide, dry chemical powder, alcohol resistant foam
- Fire fighting procedures: Wear self-contained breathing apparatus. Wear fully protective suit.

HEALTH HAZARD DATA:

Based on our knowledge, when used and handled according to specification, the product does not have harmful effects.

- Ingestion: Harmful if swallowed
- Skin contact: May cause irritation
- Eye contact: May cause irritation
- Emergency first aid procedures:
Skin: Immediately wash with water and soap. Rinse with water thoroughly.
Eyes: Flush immediately with water for at least 15 minutes. Consult doctor.
Inhaled: Supply with fresh air or oxygen. Consult doctor if discomfort persists.
Ingested: Immediately wash mouth with plenty of water. Also, call doctor immediately.

SPILL, LEAK AND DISPOSAL PROCEDURES:

If spilled or leaked, collect on absorbent. Disposal should be in accordance with ALL FEDERAL, STATE, and LOCAL ENVIRONMENTAL REGULATIONS.

PROTECTION INFORMATION:

Follow good laboratory practice when handling this product. The following precautionary measures are recommended:

- Eye protection: Laboratory safety goggles
- Hands: Chemical resistant gloves
- Skin: Laboratory protective clothing
- Respiratory: Fume hood or in areas with adequate ventilation/exhaustion.

SPECIAL PRECAUTIONS:

- Storage of sodium azide: Store at 2-8°C in well-sealed container.
- Storage of antibody containing sodium azide: Store as recommended on antibody vial, 4°C or -20°C.

This product is intended for research use only.

The above information is believed to be correct but does not purport to be all inclusive and is only intended to be used as a guide. Active Motif shall not be held responsible for any damage resulting from the use of these products or the information contained in this Material Safety Data Sheet.

Document Prepared Date: March 2005