

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.

- Stop solutions from all TransAM Kits, part #s 100347 and 100553

HAZARDOUS INGREDIENTS:

<u>Materials</u>	1N Sulphuric acid CAS# 7664-93-9
<u>Toxicity Data</u>	LD ₅₀ (Species and Route): Not applicable LD ₅₀ (Specify Species): Not applicable

The concentration of Sulphuric acid 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 DATA:

- Physical State: Liquid
- Odour and Appearance: Colorless to dark brown oily solution
- Odour threshold (ppm): Not applicable
- Boiling point: Not applicable
- Freezing point: Not applicable
- Vapor pressure: Not applicable
- Vapor density: Not applicable
- Evaporation rate: Not applicable

REACTIVITY DATA:

- Chemical Stability: Yes, stable
- Incompatible with other substances: Yes, incompatible with water and strong bases

- Reactivity: Not applicable
- Hazardous Decomposition Products: Not applicable

FIRE AND EXPLOSIONHAZARD 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: Use dry chemical or carbon dioxide. Do not use water.
- Fire fighting procedures: Wear self-contained breathing apparatus. Wear fully protective suit.

HEALTH HAZARD DATA:

Effects of acute exposure to products: Contact with skin or eyes may cause severe irritation or burns. Vapors may be irritating to eyes, nose and throat. Poison may be fatal if swallowed, causes burns.

- Effect of chronic Exposure to Product: Not applicable
- Irritancy: Not applicable
- Sensitization to Product: Not applicable
- Carcinogenicity: Not applicable
- Teratogenicity: Not applicable
- Reproductive Toxicity: Not applicable
- Mutagenicity: Not applicable
- Synergistic Products: Not applicable

PREVENTATIVEMEASURES:

- Personal Protective Equipment: This is a laboratory-use product for which no industrial protective equipment has been designed
- Gloves (specify): Latex Gloves
- Respirator (Specify): Not applicable
- Eye (Specify): Safety Glasses
- Footwear (Specify): Not applicable
- Clothing (Specify): Lab coat and apron
- Other (Specify): Not applicable
- Engineering Controls: Use with adequate ventilation
- Leak and Spill Procedure: Neutralize spill with soda ash or lime. Carefully place material into spill control container and remove from the area. Flush spill area with water.
- Waste Disposal: Dispose of in accordance with all applicable federal, state and local environmental regulations.

- Handling Procedure and Equipment: Do not get in eyes or skin or clothing. Do not breathe vapor. Wash thoroughly after handling.
- Storage Requirements: Keep in tightly closed container. Store separately and away from flammables and combustible materials.
- Special Shipping Information: Not applicable

FIRST AID MEASURES:

Specific Measures: If swallowed, **do not** induce vomiting; if conscious; give water, milk or milk of magnesia. In case of contact, immediately flush eyes or skin with plenty of water for at least 15 minutes while removing contaminated clothing and shoes. Wash clothes before re-use.

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