



MATERIAL SAFETY DATA SHEET

1 PRODUCT AND COMPANY IDENTIFICATION

Product Name: 40nmole dA(Bz) SUPPORT

Product No.: 360371; 360987; 4304078C; 4304081C; 4304079C; 4304082C; 360372; 360373; 360374; 360643; 360798; 361071; 360799; 360800; 361068; 361069; 361070; 361072; GEN759220; 4316855; 4316857; 4316859; 4316861; 4316863; 360792; 360793; 360794; 360795; 4346866

Trade Name(s): DMT-TEG POLYSTYRENE SUPPORT; HI-LOAD dA(Bz) PS SUPPORT; HI-LOAD dC(Bz) PS SUPPORT; HI-LOAD dG(iBu) PS SUPPORT; HI-LOAD dG(dmf) PS SUPPORT; HI-LOAD T PS SUPPORT; dA(Bz)SUPPORT; dC(Bz)SUPPORT; dG(iBu)SUPPORT; T SUPPORT; dG(DMF) SUPPORT; dA(Bz) 1 STEP SPRT; dC(Bz) 1 STEP SPRT; dG(iBu) 1 STEP SPRT; T 1 STEP SPRT; dG(DMF) 1 STEP SPRT; 40nMole dG (DMF) w/o OIL; 40nmole dC(Bz) w/o OIL; 40nmole 'T' w/o OIL

Manufacturer Name:

Applied Biosystems
850 Lincoln Centre Drive
Foster City, CA 94404
USA

Phone: (650) 638-5800 (USA)
www.appliedbiosystems.com

Emergency Telephone:

1-800-424-9300 (NORTH AMERICA); +1-703-527-3887 (INTERNATIONAL, REVERSE CHARGES)

Intended Use: For Research Use Only, Not for Use in Diagnostic Procedures

2 HAZARDS IDENTIFICATION

Emergency Overview

Physical State: Solid

Color: No data available.

Odor: No data available.

The toxicological properties of this product have not been thoroughly investigated. Use appropriate precautions.

Potential Health Effects

Inhalation: High concentrations: Dust may irritate the eyes and the respiratory system.

Eye Contact: Dust may irritate the eyes.

Skin Contact: No data available. However, sensitive individuals may experience slight irritation.

Ingestion: No data available. May cause discomfort if swallowed.

Target Organ(s): | Eye | Skin |

Potential Physical / Chemical Effects: Combustible solid.

OSHA Regulatory Status: This product is not hazardous according to OSHA 29CFR 1910.1200.

Environment: The product components are not classified as environmentally hazardous. However, this does not exclude the possibility that large or frequent spills can have a harmful or damaging effect on the environment.

3	COMPOSITION / INFORMATION ON INGREDIENTS
----------	-------------------------------------------------

General Information: The product contains: Reagent.

The components are not hazardous or are below required disclosure limits.

4	FIRST AID MEASURES
----------	---------------------------

Inhalation: Move injured person into fresh air and keep person calm under observation. Get medical attention if any discomfort continues.

Eye Contact: Immediately flush with plenty of water for up to 15 minutes. Remove any contact lenses and open eyes wide apart. Get medical attention if irritation develops and persists.

Skin Contact: Remove contaminated clothes and rinse skin thoroughly with water for at least 15 minutes. Get medical attention if irritation develops and persists.

Ingestion: Rinse mouth thoroughly with water and give large amounts of milk or water to people not unconscious. Get medical attention if any discomfort occurs.

5	FIRE-FIGHTING MEASURES
----------	-------------------------------

Extinguishing Media: Extinguish with foam, carbon dioxide, dry powder or water fog.

Unsuitable Extinguishing Media: Not applicable.

Special Fire Fighting Procedures: Use standard firefighting procedures and consider the hazards of other involved materials.

Unusual Fire & Explosion Hazards: Risk of dust explosion:

Hazardous Combustion Products: Carbon Dioxide, Carbon Monoxide

Protective Measures: Selection of respiratory protection for fire fighting: follow the general fire precautions indicated in the workplace. Self-contained breathing apparatus and full protective clothing must be worn in case of fire.

Flammability Class: NFPA Rating Fire = 1.

6 ACCIDENTAL RELEASE MEASURES

Personal Precautions: Avoid inhalation of dust and contact with skin and eyes. Wear suitable protective clothing. See Section 8 of the MSDS for Personal Protective Equipment.

Spill Cleanup Methods: Collect dust using a vacuum cleaner equipped with HEPA filter. For waste disposal, see section 13 of the MSDS.

Notification Procedures: Inform authorities if large amounts are involved.

7 HANDLING AND STORAGE

Handling: Keep the workplace clean. Use work methods which minimize dust production. Do not breathe dust. Avoid contact with skin and eyes. Observe good industrial hygiene practices. See Section 8 of the MSDS for Personal Protective Equipment.

Storage: Store in closed original container in a dry place. Keep away from food, drink and animal feeding stuffs.

8 EXPOSURE CONTROLS / PERSONAL PROTECTION

Exposure Limits: No exposure limits noted for ingredient(s) on ACGIH, OSHA and WEEL lists. Consult Canadian Provincial Regulations and/or Mexican Regulations on exposure limits, if applicable. However, ACGIH has limits for "nuisance dusts" which is TLV-TWA = 10 mg/m³.

Engineering Controls: Ventilate as needed to control airborne dust.

Respiratory Protection: If engineering controls do not maintain airborne concentrations below recommended exposure limits (where applicable) or to an acceptable level (in countries where exposure limits have not been established), an approved respirator must be worn. In the United States of America, if respirators are used, a program should be instituted to assure compliance with OSHA Standard 63 FR 1152, January 8, 1998. Seek advice from local supervisor.

Eye Protection: Risk of contact: Wear approved safety goggles.

Hand Protection: Wear protective gloves. Suitable gloves can be recommended by the glove supplier.

Skin Protection: Risk of contact: Wear suitable protective clothing.

Hygiene Measures: Wash hands after handling.

9 PHYSICAL AND CHEMICAL PROPERTIES

Color: No data available.

Odor: No data available.

Odor Threshold: No data available.

Physical State: Solid

pH: Not applicable

Melting Point: No data available.

Freezing Point: No data available.

Boiling Point: Not applicable.
Flash Point: No data available.
Evaporation Rate: No data available.
Flammability Limit - Upper (%): No data available.
Flammability Limit - Lower (%): No data available.
Vapor Pressure: No data available.
Vapor Density (Air=1): No data available.
Specific Gravity: No data available.
Solubility in Water: No data available.
Solubility (Other): No data available.
Partition Coefficient (n-Octanol/water): No data available.
Autoignition Temperature: No data available.
Decomposition Temperature: No data available.
Viscosity: Not applicable.

10	STABILITY AND REACTIVITY
-----------	---------------------------------

Stability: This product is stable under expected conditions of use.

Conditions to Avoid: Flame, Heat, Sparks

Incompatible Materials: Strong oxidizing agents.

Hazardous Decomposition Products:

At Elevated Temperatures:	Carbon Dioxide, Carbon Monoxide
---------------------------	---------------------------------

11	TOXICOLOGICAL INFORMATION
-----------	----------------------------------

Listed Carcinogens: None.

Product Information

Acute Toxicity: The toxicological properties of this product have not been thoroughly investigated. Use appropriate precautions.

Chronic Toxicity: No data available.

12	ECOLOGICAL INFORMATION
-----------	-------------------------------

Ecotoxicity: The product components are not classified as environmentally hazardous. However, this does not exclude the possibility that large or frequent spills can have a harmful or damaging effect on the environment.

Mobility: The product contains substances which are insoluble in water and which sediment in water systems.

Persistence and Degradability: No data available.

Bioaccumulation Potential: No data available.

13	DISPOSAL CONSIDERATIONS
-----------	--------------------------------

General Information: Dispose of waste and residues in accordance with local authority requirements.

Disposal Methods: No specific disposal method required.

Container: Since emptied containers retain product residue, follow label warnings even after container is emptied.

14	TRANSPORT INFORMATION
-----------	------------------------------

DOT Non hazardous for transport.

TDG Non hazardous for transport.

IATA Non hazardous for transport.

IMDG Non hazardous for transport.

15	REGULATORY INFORMATION
-----------	-------------------------------

Canadian Controlled Products Regulations: This product has been classified according to the hazard criteria of the Canadian Controlled Products Regulations, Section 33, and the MSDS contains all required information.

WHMIS Classification: This is not a WHMIS controlled product.

Mexican Dangerous Statement: This product is not dangerous according to Mexican regulations.

Inventory Status: US TSCA inventory Status: All the components of this substance are listed on or are exempt from the inventory. For purposes of 40 CFR 720.36, this product is for Research and Development (R&D) Use Only.

US Regulations

CERCLA Hazardous Substance List (40 CFR 302.4): Not regulated.

SARA Title III

Section 302 Extremely Hazardous Substances (40 CFR 355, Appendix A): Not regulated.

Section 311/312 (40 CFR 370):

☐ Acute (Immediate) ☐ Chronic (Delayed) ☐ Fire ☐ Reactive ☐ Pressure Generating

Section 313 Toxic Release Inventory (40 CFR 372): Not regulated.

Clean Air Act (CAA) Section 112(r) Accidental Release Prevention (40 CFR 68.130):
Not regulated.

Clean Water Act Section 311 Hazardous Substances (40 CFR 117.3): Not regulated.

Drug Enforcement Act: Not regulated.

TSCA:

TSCA Section 4(a) Final Test Rules & Testing Consent Orders: Not regulated.

TSCA Section 5(a)(2) Final Significant New Use Rules (SNURs) (40CFR 721, Subpt. E): Not regulated.

TSCA Section 5(e) PMN-Substance Consent Orders: Not regulated.

TSCA Section 12(b) Export Notification (40 CFR 707, Subpt. D): Not regulated.

State Regulations

California Safe Drinking Water and Toxic Enforcement Act of 1986 (Proposition 65):
Not regulated.

Massachusetts Right-To-Know List: Not regulated.

Michigan Critical Materials List (Michigan Natural Resources and Environmental Protection Act (Act. 451 of 1994)): Not regulated.

Minnesota Hazardous Substances List: Not regulated.

New Jersey Right-To-Know List: Not regulated.

Pennsylvania Right-To-Know List: Not regulated.

Rhode Island Right-To-Know List: Not regulated.

16	OTHER INFORMATION
-----------	--------------------------

HAZARD RATINGS

	Health Hazard	Fire Hazard	Instability	Special Hazard
NFPA	1	1	0	--

Hazard rating: 0 - Minimal; 1 - Slight; 2 - Moderate; 3 - Serious; 4 - Severe

NFPA Label colored diamond code: Blue - Health; Red - Flammability; Yellow - Instability; White - Special Hazards

	Health Hazard	Flammability	Physical Hazard	Personal Protection
HMIS	1	1	0	B

Hazard rating: 0 - Minimal; 1 - Slight; 2 - Moderate; 3 - Serious; 4 - Severe

Personal Protection codes: B - Safety Glasses, Gloves

HMIS Label colored bar code: Blue - Health; Red - Flammability; Orange - Physical Hazards; White - Special

General Information: To the best of our knowledge, the chemical, physical, and toxicological properties of this material have not been thoroughly investigated. The potential hazards described in this MSDS are based on the structure activity relationship of structurally similar chemicals, or similar chemicals belonging to the same chemical class for which data is available.

Disclaimer: This MSDS has been prepared using information from sources considered technically reliable. It should not be relied upon as a product specification. The company makes no warranty of any kind, express or implied, concerning the safe use of this material in your process or in combination with other chemical substances. Every user of the product is responsible for evaluating its own processes and conditions of use and selecting appropriate protective measures for employee exposure.