Safety Data Sheet (SDS)

Revision Number: 3.0		Last updated March 17, 2021	
1. Product and Company Ider	ntification_		
Product Name:	5(6) - TRIT	5(6) - TRITC	
	[Tetramethy	rrhodamine - 5 - (and - 6) - isothiocyanate	
Manufacturer/Supplier:	AnaSpec, In	ic.	
	www.anasp	ec.com	
	34801 Cam	pus Drive	
	Fremont, CA 94555		
	Tel: 510-791-9560		
	Fax: 510-791-9572		
	Email: servi	ce@anaspec.com	
	Kaneka Eur	ogentec SA,	
	Rue du Bois	s Saint Jean 5 4102 Seraing Belgium	
	Tel. +32-4-3	3727400	
	Fax. +32-4-	3727500	
	E-mail info	@eurogentec.com	
	Kaneka Eur	ogentec Helpdesk	
	Tel. +32-4-3	•	
Catalog Number	AS-81150		

2. Hazards Identification

Emergency Overview: We do recommend handling all chemicals with caution. Use proper protective equipment when handling chemicals. To our knowledge, the hazards of this material have not been thoroughly investigated.

GHS Hazard Classification:

Skin irritation (Category 2) Eye irritation (Category 2A)

Respiratory sensitization (Category 1)

Skin sensitization (Category 1)

Specific target organ toxicity - single exposure (Category 3)

GHS Signal Words: DANGER

GHS Hazard Symbol/Pictogam:



GHS Hazard Statements:

H315 Causes skin irritation

H317 May cause an allergic skin reaction.

H319 Causes serious eye irritation

H334 May cause allergy or asthma symptoms or breathing difficulties if inhaled

H335 May cause respiratory irritation

GHS Precautionary Statements:

P261 Avoid breathing dust/ fume/ gas/ mist/ vapours/ spray

P280 Wear protective gloves

P305+P351+P338 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact

lenses, if present and easy to do. Continue rinsing

P342+P311 If experiencing respiratory symptoms: Call a POISON CENTER or doctor/

physician

Potential Health Effects for:

Inhalation: If inhaled, remove to fresh air. If not breathing, give artificial respiration. If breathing is difficult, give

oxygen. Good hygiene practice requires that exposure be kept to a minimum and that suitable control

measures be used in an occupational setting.

Ingestion: If swallowed, wash out mouth with water provided person is conscious. Call a physician.

Skin: In case of contact, immediately wash skin with soap and copious amount of water.

Eyes: In case of contact, immediately flush eyes with copious amounts of water for at least 15 minutes.

Chronic Exposures: No information available. We recommend limiting prolonged exposure.

Target Organs: No information available

HMIS Classification

Health hazard: 1
Flammability: 0
Reactivity Hazard: 0

NFPA Rating

Health hazard: 1

Fire: 0

Reactivity Hazard: 0

3. Composition / Information on Ingredients

Chemical Name: 5(6)-TRITC; [Tetramethylrhodamine –5-(and 6)-isothiocyanate];

xanthylium, 9-(2-carboxy isothiocyantophenyl)-3,6-bis-(dimethylamino)-inner salt

Molecular Formula: C25H21N3O3S

Molecular Weight: 443.5 CAS –No: 95197-95-8

Ingredients/C	Components:		
4. First Aid	Measures		
Inhalation:			
		ent to blow nose to ensure clear passage of breathing.	
	If irritation or discomfort persists seek medical attention.		
Ingestion: If swallowed do NOT inc If vomiting occurs, lean p maintain open airway and Observe the patient caref			
		occurs, lean patient forward or place on left side (head-down position, if possible) to	
		nse out mouth, then provide liquid slowly and as much as casualty can comfortably	
	drink.	,,,	
	Seek medical ad	lvice	
Skin:	If skin or hair c		
		nair with running water (and soap if available).	
E		tention in event of irritation.	
Eyes:		omes in contact with the eyes:	
		nediately with fresh running water. ete irrigation of the eye by keeping eyelids apart and away from eye and moving the	
		sionally lifting the upper and lower lids.	
		or recurs seek medical attention.	
	-		
5. Fire Fight	ting Measures		
Extinguishing	media:	Water spray or fog.	
		Alcohol resistant foam.	
		Dry chemical powder.	
		BCF (where regulations permit). Carbon dioxide	
Special firefic	ghting procedures:	Alert Emergency Responders and tell them location and nature of	
<i>Бресіш</i> зігезія	gning procedures.	hazard.	
		Wear breathing apparatus plus protective gloves.	
		Prevent, by any means available, spillage from entering drains or water	
		course.	
		Use water delivered as a fine spray to control fire and cool adjacent	
		area.	
		DO NOT approach containers suspected to be hot.	
		Cool fire exposed containers with water spray from a protected	
		location.	
		If safe to do so, remove containers from path of fire. Equipment should be thoroughly decontaminated after us	
Unusual fire	and explosions haz		
Onusuai jire t	and explosions haz	areas. Elines toxic funcional fire conditions	
6. Accidenta	ıl Release Measuro	<u>es</u>	
Spill response	 2	Remove all ignition sources.	
F 122p 0.050		Clean up all spills immediately.	
		avoid contact with skin and eyes.	
		Control personal contact by using protective equipment.	
		Jse dry clean up procedures and avoid generating dust.	
		Place in a suitable, labeled container for waste disposal	
Containment	Avoid all personal contact, including inhalation.		

Wear protective clothing when risk of exposure occurs.

Use in a well-ventilated area.

DO NOT enter confined spaces until atmosphere has been checked.

DO NOT allow material to contact humans, exposed food or food utensils.

Avoid contact with incompatible materials.

When handling, DO NOT eat, drink or smoke.

Keep containers securely sealed when not in use.

Avoid physical damage to containers.

Always wash hands with soap and water after handling.

Use good occupational work practice.

Empty containers may contain residual dust which has the potential to accumulate following settling. Such dusts may explode in the presence of an appropriate ignition source.

Do NOT cut, drill, grind or weld such containers

PPE

Use personal protective equipment

7. Handling and Storage

Store at 4 °C desiccated and protected from light. Store away from oxidizing agent.

8. Exposure Controls / Personal Protection

			. 7
Hn	ain.	ooring	controls
LIL	zuw	ceinig	Commons

Local exhaust ventilation is required where solids are handled as powders or crystals; even when particulates are relatively large, a certain proportion will be powdered by mutual friction.

Exhaust ventilation should be designed to prevent accumulation and re-circulation of particulates in the workplace.

If in spite of local exhaust an adverse concentration of the substance in air could occur, respiratory protection should be considered. Such protection might consist of:

(a): particle dust respirators, if necessary, combined with an absorption cartridge;

(b): filter respirators with absorption cartridge or canister of the right type:

(c): fresh-air hoods or masks

Build-up of electrostatic charge on the dust particle, may be prevented by bonding and grounding.

Powder handling equipment such as dust collectors, dryers and mills may require additional protection measures such as explosion venting.

Air contaminants generated in the workplace possess varying "escape" velocities which, in turn, determine the "capture velocities" of fresh circulating air required to efficiently remove the contaminant.

PPE

Respiratory protection

Where risk assessment shows air-purifying respirators are appropriate use a full-face respirator with multi-purpose combination (US) or type ABEK (EN 14387) respirator cartridges as a backup to engineering controls. If the respirator is the sole means of protection, use a full-face supplied air respirator. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

Hand protection

Handle with gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.

Eye protection

AnaSpec Inc.

Toll-Free: 800-452-5530 • Tel: 510-791-9560 • Fax: 510-791-9573

Face shield and safety glasses Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU). Skin and body protection

Complete suit protecting against chemicals, Flame retardant antistatic protective clothing, The type of protective equipment must be selected according to the concentration and amount of the dangerous substance at the specific workplace.

Hygiene measures

Avoid contact with skin, eyes and clothing. Wash hands before breaks and immediately after handling the product

9. Physical and Chemical Properties

Physical State	Solution
Odour	Not available
Solubility in Water	Not available
Specific Gravity	Not available
рН	Not available
Boiling Point	N/A
Melting Point	Not available
Flash Point	N/A
Vapor Pressure:	N/A
Vapor Density:	N/A

10. Stability and Reactivity

Thermal Decomposition	No data available
Dangerous Products of Decomposition	No data available
Dangerous Reactions	COx, NOx when burned

Keep container tightly closed in a dry well-ventilated place. Store in 4°C refrigerator.

11. Toxicological Information

RTECS Number	N/A
Toxicity	No information available.
Health Hazards	Although ingestion is not thought to produce harmful effects, the material may still be damaging to the health of the individual following ingestion, especially where pre-existing organ (e.g. liver, kidney) damage is evident. In an occupational setting however, ingestion of insignificant quantities is not thought to be cause for concern.
Potential Hazards	Not available
Carcinogenicity:	No significant acute toxicological data identified
OSHA Permissible Exposure Limit(PEL) Data	N/A
ACGIH Threshold Limit Values (TLV)	N/A
Panyaduatina Tariaitu	No information available

Reproductive Toxicity: No information available

12. Ecological Information

No information available.

13. Disposal Considerations

All waste must be handled in accordance with local, state and federal regulations. Legislation addressing waste disposal requirements may differ by country, state and/ or territory. Each user must refer to laws operating in their area. In some areas, certain wastes must be tracked.

14. Transport Information

Hazard Class	N/A
Identification Number	N/A
Packing Group	N/A
Proper Shipping Name (DOT)	N/A

15. Regulatory Information

California Proposition 65: N/A

US TSCA (Toxic Substance Control Act): N/A

US CERCLA (Comprehensive Environmental Response, Compensation, and Liability Act: N/A

US SARA Title III (Superfund Amendments and Reauthorization Act: N/A

US Other: N/A

EC EINICS (European Inventory of Existing Commercial Chemical Substances) Number: N/A

EC Risk Statements: N/A

Other Country Regulations: N/A

16. Other Information

It is not intended for food, drug, household, agricultural or cosmetic use. A technically qualified individual experienced in handling potentially hazardous chemicals must supervise its use. The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. Users should make independent decisions regarding completeness of the information based on all sources available. AnaSpec shall not be held liable for any damage resulting from handling or from contact with the above product.