Safety Data Sheet (SDS)

Revision Number: 3.0		Last updated	9 April 2020
1. Product and Company Identification	<u>on</u>		
Product Name:	Spike protein (S1/S2) SARS - CoV - 2 substrate QXL® 520 - TNS PRR ARS VAS K(5-FAM) - NH2		
Manufacturer/Supplier:	AnaSpec, Inc. www.anaspec.com 34801 Campus Drive Fremont, CA 94555 Tel: 510-791-9560 Fax: 510-791-9572 Email: service@anaspec.com Kaneka Eurogentec SA, Rue du Bois Saint Jean 5 4102 Seraing Belgium Tel. +32-4-3727400 Fax. +32-4-3727500 E-mail info@eurogentec.com		
	Kaneka Eur Tel. +32-4-3	ogentec Helpdesk 3727665	
Catalog Number	AS-65600		
Relevant identified uses of the substance/preparation and uses advised against	For laborate	ry use only.	
Emergency information		act the regional Eurogentec representation and Eurogentec S.A. directly (from	•
2. Hazards Identification			
		dling all chemicals with caution. Use proper of this materials. To our knowledge, the hazards of this materials.	
	-	ance according to the GHS angerous substance according to the GHS	
GHS Signal Words: None			

GHS Hazard Statements: None

GHS Precautionary Statements: None

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: 0

Chronic Health Hazard: 0

Flammability: 0
Physical hazards: 0

NFPA Rating

Health hazard: 0

Fire: 0

Reactivity Hazard: 0

3. Composition

Ingredients/Components:

Chemical Name: Spike protein (S1/S2) SARS - CoV - 2 substrate

QXL® 520 - TNS PRR ARS VAS K(5-FAM) - NH2

Molecular formula: NA Molecular weight: 2278.4

CAS-No NA EC-No NA

4. First Aid Measures

Inhalation:	If dust is inhaled, remove from contaminated area.
	Encourage patient to blow nose to ensure clear passage of breathing.
	If irritation or discomfort persists seek medical attention.
Ingestion:	If swallowed do NOT induce vomiting.
	If vomiting occurs, lean patient forward or place on left side (head-down position, if possible) to maintain open airway and prevent aspiration.
	Observe the patient carefully.
	Give water to rinse out mouth, then provide liquid slowly and as much as casualty can comfortably
	drink.
	Seek medical advice.
Skin:	If skin or hair contact occurs:
	Flush skin and hair with running water (and soap if available).
	Seek medical attention in event of irritation.
Eyes:	If this product comes in contact with the eyes:
	Wash out immediately with fresh running water.
	Ensure complete irrigation of the eye by keeping eyelids apart and away from eye and moving the
	eyelids by occasionally lifting the upper and lower lids.
	If pain persists or recurs seek medical attention.

T		XXI.	
Extinguishing media:		Water spray or fog.	
		Alcohol resistant foam.	
		Dry chemical powder. BCF (where regulations permit).	
		Carbon dioxide	
		curson dismue	
Special firefighting procedures:		Alert Emergency Responders and tell them location and nature of 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 use.	
Unusual fire and explosions hazards:		Emits toxic fumes under fire conditions	
6. Accidental Release	<u>Measures</u>		
Spill response		ll ignition sources.	
		all spills immediately.	
		ntact with skin and eyes.	
		ersonal contact by using protective equipment.	
		ean up procedures and avoid generating dust. suitable, labeled container for waste disposal	
Containment		personal contact, including inhalation.	
	Wear prot	ective clothing when risk of exposure occurs.	
		vell-ventilated area.	
		enter confined spaces until atmosphere has been checked.	
		allow material to contact humans, exposed food or food utensils.	
When hat Keep con		ntact with incompatible materials. dling, DO NOT eat, drink or smoke.	
		ainers securely sealed when not in use.	
		id physical damage to containers.	
		Always wash hands with soap and water after handling.	
		occupational work practice.	
	Empty con	ntainers may contain residual dust which has the potential to accumulate	
	_	settling. Such dusts may explode in the presence of an appropriate ignition	
	source. Do NOT o	cut, drill, grind or weld such containers	
PPE	Use perso	Use personal protective equipment	

8 Exposure Controls	/ Personal Protection		
8. Exposure Controls Engineering controls	Local exhaust ventilate even when particulate mutual friction. Exhaust ventilation separticulates in the word of the separticulates (a): particle dust respirators (b): filter respirators (c): fresh-air hoods of the separticulates of the separticulates (c): fresh-air hoods (c): fresh-	haust an adverse concentration of the substance in air could occur, in should be considered. Such protection might consist of: birators, if necessary, combined with an absorption cartridge; with absorption cartridge or canister of the right type; or masks attic charge on the dust particle, may be prevented by bonding and aipment such as dust collectors, dryers and mills may require measures such as explosion venting. The relation of the substance in air could occur, and in the workplace possess varying "escape" velocities which, as "capture velocities" of fresh circulating air required to efficiently	
	remove the contamin		
PPE	Use personal protecti		
9. Physical and Chemi	cal Properties		
Physical State	Solid		
Odour	Not available		
Solubility in Water	Not available		
Specific Gravity	Not available		
pН	Not available		
Boiling Point	Not available		
Melting Point	Not available		
Flash Point	N/A		
Vapor Pressure:	N/A		
Vapor Density:	N/A		
10. Stability and Read	etivit <u>y</u>		
Thermal Decomposition	i	No data available	
*		No data available	
		COx, NOx when burned	
Keep container tightly of	closed in a dry well-ven	ntilated place. Store in -20 °C, dry refrigerator.	
11. Toxicological Info	<u>rmation</u>		
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

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.