

Material Safety Data Sheet

V1, Reviewed on: 13/10/2017
Printing date: 13/10/2017

1. Identification of the Product and of the company/undertaking

1.1 Product identifiers

- Product type: RNA Library Preparation Kit
- Product name: CATS RNA-seq Kit v2 x12, x24, x96
- Product number: C05010045, C05010041, C05010049
- REACH No.: Not available

1.2 Relevant identified uses of the substance or mixture and used-s advised against
identified uses: RUO laboratory chemicals

1.3. Details of the supplier of the safety data sheet

- Manufacturer/Supplier: Diagenode S.A.
Liège Science Park
Rue du Bois St-Jean, 3
4102 Seraing (Ougrée)
Belgium
- Tel: +32-4-364-20-50
- Fax: +32-4-364-20-51
- email: www.diagenode.com

1.4. Information in case of emergency:

Please contact the regional Diagenode representation in your country or Diagenode S.A. directly. For a complete listing of Diagenode's international distributors visit: <http://www.diagenode.com/pages/distributors.html>

2. Hazards identification

2.1. Classification of the substance or mixture

- Acute toxicity, category 4, H302
- Serious eye damage, category 1, H318
- Eye irritation, category 2, H319
- Skin irritation, category 2, H315
- Chronic aquatic toxicity (Category 2), H411
- Classification system: The classification is according to the latest editions of the EU-lists, and extended by company and literature data.

2.2. Label elements

pictogram



signal word

warning

hazard statements

- H302: harmful if swallowed
- H315: causes skin irritation
- H318: causes serious eye damage
- H319: causes serious eye irritation
- H335: may cause respiratory irritation
- H411: toxic to aquatic life with long-lasting effects

precautionary statements

- P261: Avoid breathing dust/fume/gas/mist/vapours/spray
- P280: Wear protective gloves/protective clothing/eye protection/face protection
- P301 + P312 + P330 IF SWALLOWED: Call a POISON CENTER/doctor if you feel unwell. Rinse mouth.
- P305 + P351 + P338 + P310 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Immediately call a POISON CENTER/doctor

2.3. Other hazards

None

3. Composition/information on ingredients

3.1. Substances

Not applicable

3.2. Mixtures

This product is composed of buffers and synthetic DNA oligonucleotides in aqueous buffer solutions. It contains the following potentially hazardous chemicals:

Chemical name	CAS No.	EC No.	Classification	Concentration
Triton X-100	9002-93-1	618-344-0	Acute Tox. 4, H302; Skin Irrit. 2, H315; Eye Dam. 1, H318; Aquatic Chronic 2, H411	≤0.2%
Dithiothreitol (DTT)	3483-12-3	222-468-7	Acute Tox. 4, H302; Skin Irrit. 2, H315; Eye Irrit. 2, H319; STOT SE 3, H335	≤0.2%

4. First-aid measures

4.1. Description of first aid measures

- **After inhalation**

Provide fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen. Get medical attention.

- **After skin contact**

Flush contaminated areas with water for at least 15 minutes. If irritation persists, get medical attention.

- **After eye contact**

Rinse eyes with plenty of water for at least 15 minutes. If irritation persists, get medical attention.

- **After swallowing**

Rinse mouth with water. Do NOT induce vomiting unless directed to do so by medical personnel. Never give anything by mouth to an unconscious person. Get medical attention.

4.2. Most important symptoms and effects, both acute and delayed

See section 2 and 11

4.3. Indication of any immediate medical attention and special treatment needed

No data available

5. Fire-fighting measures

5.1. Extinguishing media

Suitable extinguishing agents: Water, CO₂, powder.

5.2. Special hazards arising from the substance or mixture

No data available

5.3. Advice for fire-fighters

No data available

6. Accidental release measures

6.1. Personal precautions, protective equipment and emergency procedures

Use personal protective equipment. Avoid breathing vapours, mist or gas. Ensure adequate ventilation.

6.2. Environmental precautions

Prevent further leakage or spillage. Do not allow product to reach sewage system or any water course. Discharge into the environment must be avoided.

6.3. Methods and material for containment and cleaning up

Absorb with inert absorbent material and dispose of as hazardous waste. Keep in suitable, closed containers for disposal.

6.4. Reference to other sections

See section 8 and 13

7. Handling and storage

7.1. Precautions for safe handling

Avoid inhalation and contact with eyes and skin.

7.2. Conditions for safe storage, including any incompatibilities

Keep container tightly closed in a dry and well ventilated place. Containers which are opened should be carefully resealed and kept upright to prevent leakage.

7.3. Specific end use

See section 1.2.

8. Exposure controls/personal protection

8.1. Control parameters

No data available

8.2. Exposure controls

- **Engineering Controls:** Safety shower and eye bath. Use only in a chemical fume hood.
- **Personal Protection Equipment:** Lab coat. Compatible chemical-resistant gloves. Safety glasses. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).
- **General Hygiene Measures:** Wash contaminated clothing before reuse. Wash thoroughly after handling.

Exposure Limits:	Country	Source	Type	Value
	USA	ACGIH	TWA	Not Available
	USA	OSHA	PEL	Not Available

9. Physical and chemical properties

9.1. Information on basic physical and chemical properties

Physical State And Appearance: Liquid, clear

Chemical Formula: Not available

Boiling Point: Not available

Melting Point: Not available

Freezing Point: Not available

Flash Point: Not available

Auto Igniting: Not available

Vapor Pressure: Not available

Vapor Density: Not available

SG/Density: Not available

Flammability: Not available

Autoignition Temp: Not available

Solubility: Not available

9.2. Other information

No data available

10. Stability and reactivity

10.1. Reactivity

No data available

10.2. Chemical stability

Stable until the provided expiry date when stored under the recommended conditions

10.3. Possibility of hazardous reactions

No data available

10.4. Conditions to avoid

No data available

10.5. Incompatible materials

Avoid strong oxidizing agents.

10.6. Hazardous decomposition products

- Thermal decomposition: May form carbon dioxide and carbon monoxide.
- Others: No data available.

11. Toxicological information

11.1. Information on toxicological effects

- RTECS No. Not available
- Potential Health Effects:
 - Inhalation May be harmful if inhaled. May cause respiratory tract irritation.
 - Ingestion May be harmful if swallowed.
 - Skin May be harmful if absorbed through skin. May cause skin irritation.
 - Eyes May cause eye irritation.
- Signs And Symptoms Of Exposure:

To the best of our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated. Use appropriate procedures and precautions to prevent or minimize exposure.
- Acute Toxicity
 - Oral LD50: No data available
 - Inhalation LC50: No data available
 - Dermal LD50: No data available

12. Ecological information

12.1. Toxicity

No data available

12.2. Persistence and degradability

No data available

12.3. Bioaccumulative potential

No data available

12.4. Mobility in soil

No data available

12.5. Results of PBT and vPvB assessment

No data available

12.6. Other adverse effects

No data available

13. Disposal considerations

13.1. Waste treatment methods

- Product: Dispose as special waste in compliance with official regulations.
- Uncleaned packaging: Dispose in compliance with official regulations.

14. Transport information

14.1. UN number

Not applicable

14.2. UN proper shipping name

Not applicable

14.3. Transport hazard class(es)

ADR: Not classified as dangerous in the meaning of transport regulations.

IMDG: Not classified as dangerous in the meaning of transport regulations.

IATA: Not classified as dangerous in the meaning of transport regulations.

14.4. Packing group

Not applicable

14.5. Environmental hazards

None

14.6. Special precautions for user

No data available

14.7. Transport in bulk according to Annex II of MARPOL 73/78 and the IBC Code

Not applicable

15. Regulatory information

15.1. Safety, health and environmental regulations/legislation specific for the substance or mixture

This safety datasheet complies with the requirements of Regulation (EC) No. 1907/2006

15.2. Chemical Safety Assessment

For this product a chemical safety assessment was not carried out

16. Other information

This information is based on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship. The information in this document is intended only as a guide to the appropriate precautionary handling of this material by a person trained in, or supervised by a person trained in chemical handling. The user is responsible for determining the precautions and danger of this chemical material in his application. This material may react with other substances and since potential users are so varied, Diagenode S.A., cannot warn of all potential dangers. Therefore Diagenode S.A. will not assume responsibility and expressly disclaims liability for loss, damage, or expense arising from the use, handling, storage, or disposal of this product. Diagenode S.A. urges each user to carefully read the MSDS prior to using the product and become aware to the adverse health effects associated with this product. While the information provided in the MSDS is believed to be correct, the information is provided without any warranty expressed or implied, regarding its corrective uses.