

**Section 1. Identification**

**Product identifier**

**Product Identity** Armored RNA®, Armored RNA® Quant™, Custom Armored RNA®, Custom Armored RNA® Quant™, Armored DNA Quant™, Custom Armored DNA Quant™

**Other means of identification** 42000, 42004, 42006, 42008, 42010, 42012, 42024, 42030, 42032, 42040, 42041, 42042, 42044, 42045, 42046, 42047, 42050, 42090, 42091, 42101, 42102, 42103, 49610, 49650, 49659, 49705, 52000, 52003, 52030, 52031, 52036

**Relevant identified uses of the substance or mixture and uses advised against**

Lab reagents

**Details of the supplier of the safety data sheet**

**Company Name** Asuragen, Inc.  
2150 Woodward Street  
Austin, Texas 78744  
+1 512 681 5200

**Emergency  
24 hour Emergency  
Telephone No.** Asuragen, Inc.  
+1 512 681 5200 US, Toll-free Tel: 1-877-777-1874 (Normal Business Hours)

**Customer Service:**

## Section 2. Hazard(s) identification

Classification of the substance or mixture under US OSHA's Hazard Communication Standard (1910.1200) revised 2024 and Canadian Hazardous Products Regulations (SOR/2015-17) (GHS revision 7)

### Label elements

#### [Prevention]

No GHS prevention statements

#### [Response]

No GHS response statements

#### [Storage]

No GHS storage statements

#### [Disposal]

No GHS disposal statements

### Other hazards

This product contains no PBT/vPvB/vPvM chemicals.

This product contains no endocrine disrupting chemicals.

Does not contain component(s) that meet a 'definition' of per & poly fluoroalkyl substance (PFAS) per the Organisation for Economic Co-operation and Development (OECD) list of Per- and Polyfluoroalkyl Substances (PFASs).

## Section 3. Composition/information on ingredients

There are no ingredients in this product which are classified as hazardous, and/or no hazardous ingredients above the GHS cut off percentage.

#### Section 4. First aid measures

##### Description of first aid measures

- General** In all cases of doubt, or when symptoms persist, seek medical attention. Never give anything by mouth to an unconscious person.
- Inhalation** Remove to fresh air, keep patient warm and at rest. If breathing is irregular or stopped, give artificial respiration. If unconscious, place in the recovery position and obtain immediate medical attention. Give nothing by mouth.
- Eyes** Irrigate copiously with clean water for at least 15 minutes, holding the eyelids apart and seek medical attention.
- Skin** Remove contaminated clothing. Wash skin thoroughly with soap and water or use a recognized skin cleanser.
- Ingestion** If swallowed obtain immediate medical attention. Keep at rest. Do NOT induce vomiting.

##### Most important symptoms and effects, both acute and delayed

- Overview** No specific symptom data available.  
No chronic toxicity or long term toxicity information available. Treat symptomatically. See section 2 for further details.

#### Section 5. Fire-fighting measures

##### Extinguishing media

Recommended extinguishing media; alcohol resistant foam, CO<sub>2</sub>, powder, water spray.  
Unsuitable extinguishing media: Do not use; water jet.

**Special hazards arising from the substance or mixture**

Hazardous decomposition: No hazardous decomposition data available.

**Advice for fire-fighters**

As with all fires, wear positive pressure, self-contained breathing apparatus, (SCBA) with a full face piece and protective clothing. Persons without respiratory protection should leave area. Wear SCBA during clean-up immediately after fire. No smoking.

**ERG Guide No.**

----

**Section 6. Accidental release measures**

**Personal precautions, protective equipment and emergency procedures**

Put on appropriate personal protective equipment (see section 8).  
Use good personal hygiene practices. Wash hands before eating, drinking, smoking or using toilet. Promptly remove soiled clothing and wash thoroughly before reuse.

**Environmental precautions**

Do not allow spills to enter drains or waterways.

**Methods and material for containment and cleaning up**

Ventilate the area and avoid breathing vapors. Take the personal protective measures listed in section 8.  
Contain and absorb spillage with non-combustible materials e.g. sand, earth, and vermiculite. Place in closed containers outside buildings and dispose of according to the Waste Regulations.

**Section 7. Handling and storage**

**Precautions for safe handling**

Handle containers carefully to prevent damage and spillage.  
See section 2 for further details. - [Prevention]

**Conditions for safe storage, including any incompatibilities**

Incompatible materials: No available information  
See section 2 for further details. - [Storage]

**Specific end use(s)**

No available information

**Section 8. Exposure controls / personal protection**

There are no ingredients in this product which are classified as hazardous, and/or no hazardous ingredients above the GHS cut off percentage.

**Exposure  
controls**

<b>Respiratory</b>	If workers are exposed to concentrations above the exposure limit they must use the appropriate, certified respirators.
<b>Eyes</b>	Protective safety glasses recommended
<b>Skin</b>	Avoid skin contact. Protective gloves recommended.
<b>Engineering Controls</b>	Provide adequate ventilation. Where reasonably practicable this should be achieved by the use of local exhaust ventilation and good general extraction. If these are not sufficient to maintain concentrations of particulates and any vapor below occupational exposure limits suitable respiratory protection must be worn.
<b>Other Work Practices</b>	Use good personal hygiene practices. Wash hands before eating, drinking, smoking or using toilet. Promptly remove soiled clothing and wash thoroughly before reuse.

See section 2 for further details.

## Section 9. Physical and chemical properties

### Information on basic physical and chemical properties

<b>Physical State</b>	Liquid
<b>Color</b>	Clear
<b>Odor</b>	No available information
<b>Odor threshold</b>	No available information
<b>Melting point / freezing point</b>	No available information
<b>Initial boiling point and boiling range</b>	No available information
<b>Flammability (solid, gas)</b>	Not Applicable
<b>Upper/lower flammability or explosive limits</b>	<b>Lower Explosive Limit:</b> No available information <b>Upper Explosive Limit:</b> No available information
<b>Flash Point</b>	°F °C, Test method: (Open/Close cup)
<b>Auto-ignition temperature</b>	No available information
<b>Decomposition temperature</b>	No available information
<b>pH</b>	No available information
<b>Viscosity (cSt)</b>	No available information
<b>Solubility in Water</b>	No available information
<b>Partition coefficient n-octanol/water (Log Kow)</b>	No available information
<b>Vapor pressure (Pa)</b>	No available information
<b>Relative Density</b>	No available information
<b>Vapor Density</b>	No available information
<b>Evaporation rate (Ether = 1)</b>	No available information
<b>Oxidising properties</b>	No available information
<b>Explosive properties</b>	No available information

### Other information

No other relevant information.

## Section 10. Stability and reactivity

### Reactivity

Hazardous Polymerization will not occur.

### Chemical stability

Stable under normal circumstances.

### Possibility of hazardous reactions

No available information

### Conditions to avoid

Avoid high temperatures and contact with incompatible material

### Incompatible materials

No available information

### Hazardous decomposition products

No hazardous decomposition data available.

## Section 11. Toxicological information

### Acute toxicity

There are no ingredients in this product which are classified as hazardous, and/or no hazardous ingredients above the GHS cut off percentage.

Classification	Category	Hazard Description
Acute toxicity (oral)	---	Not Applicable
Acute toxicity (dermal)	---	Not Applicable
Acute toxicity (inhalation)	---	Not Applicable
Skin corrosion/irritation	---	Not Applicable
Serious eye damage/irritation	---	Not Applicable
Respiratory sensitization	---	Not Applicable
Skin sensitization	---	Not Applicable
Germ cell mutagenicity	---	Not Applicable

Carcinogenicity	---	Not Applicable
Reproductive toxicity	---	Not Applicable
STOT-single exposure	---	Not Applicable
STOT-repeated exposure	---	Not Applicable
Aspiration hazard	---	Not Applicable

**Possible routes of entry:**No available information

**Symptoms and effects, both acute and delayed:**

No specific symptom data available.

No chronic toxicity or long term toxicity information available. Treat symptomatically.

**Section 12. Ecological information**

**Toxicity**

No additional information provided for this product. See Section 3 for chemical specific data.

**Aquatic Ecotoxicity**

There are no ingredients in this product which are classified as hazardous, and/or no hazardous ingredients above the GHS cut off percentage.

**Persistence and degradability**

There is no data available on the preparation itself.

**Bioaccumulative potential**

No available information

**Mobility in soil**

No available information

**Results of PBT and vPvB assessment**

This product contains no PBT/vPvB/vPvM chemicals.

**Other adverse effects**

No available information



### Section 13. Disposal considerations

#### Waste treatment methods

Observe all federal, provincial and local regulations when disposing of this substance.

### Section 14. Transport information

Classification Method: Classified as per Part 2, Sections 2.1-2.8 of the Transportation of Dangerous Goods Regulations.

	<b>DOT (Domestic Surface Transportation)</b>	<b>TDG (Domestic Surface Transportation)</b>
<b>UN number</b>	Not Regulated	Not Regulated
<b>UN proper shipping name</b>	Not Regulated	Not Regulated
<b>Transport hazard class(es)</b>	<b>Class:</b> Not Applicable	<b>Class:</b> Not Applicable <b>Sub Class:</b> Not Applicable
<b>Packing group</b>	Not Applicable	Not Applicable
	<b>IMO / IMDG (Ocean Transportation)</b>	<b>ICAO/IATA</b>
<b>UN number</b>	Not Regulated	Not Regulated
<b>UN proper shipping name</b>	Not Regulated	Not Regulated
<b>Transport hazard class(es)</b>	<b>Class:</b> Not Applicable <b>Sub Class:</b> Not Applicable	<b>Class:</b> Not Applicable <b>Sub Class:</b> Not Applicable
<b>Packing group</b>	Not Applicable	Not Applicable

#### Environmental hazards

IMDG Marine Pollutant: No;

**Special precautions for user**

No available information

**Section 15. Regulatory information**

**Regulatory  
Overview**

The regulatory data in Section 15 is not intended to be all-inclusive, only selected regulations are represented.

This product has been classified in accordance with US OSHA's Hazard Communication Standard (1910.1200) revised 2024 and Canadian Hazardous Products Regulations (SOR/2015-17 amended 2022-12-15) (GHS revision 7) and the SDS contains all of the information required by those regulations.

**Toxic Substance Control Act (TSCA)**

2-methyl-3-2H-isothiazolone (PMN; SP)

5-Chloro-2-methyl-2H-isothiazol-3-one (PMN; SP)

Gelatin (UVCB XU)

Magnesium nitrate

Sodium chloride

Tromethamine

Water

**EPCRA 302 Extremely Hazardous:**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**EPCRA 313 Toxic Chemicals:**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Canadian Domestic Substance List (DSL):**

2-methyl-3-2H-isothiazolone

5-Chloro-2-methyl-2H-isothiazol-3-one

Deoxyribonucleic acids

Gelatin

Magnesium nitrate

Sodium chloride

Tromethamine

Water

**Canadian Non-Domestic Substance List (NDSL):**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Proposition 65 - Carcinogens (>0.0%):**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Proposition 65 - Developmental Toxins (>0.0%):**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Proposition 65 - Female Repro Toxins (>0.0%):**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Proposition 65 - Male Repro Toxins (>0.0%):**

To the best of our knowledge, there are no chemicals at levels which require reporting under this statute.

**Proposition 65 Label Warning:**

This product contains no chemicals known to the State of California to cause cancer and birth defects or other reproductive harm.

<b>Section 16. Other information</b>
--------------------------------------

**Revision Date**

08/05/2025

The information and recommendations contained herein are based upon data believed to be correct. However, no guarantee or warranty of any kind, expressed or implied, is made with respect to the information contained herein. We accept no responsibility and disclaim all liability for any harmful effects which may be caused by exposure to our products. Customers/users of this product must comply with all applicable health and safety laws, regulations, and orders.

The full text of the phrases appearing in section 3 is:

Not Applicable

Disclaimer: The information presented herein is supplied as a guide to those who handle or use this product. Safe work practices must be employed when working with any materials. It is important that the end user makes a determination regarding the adequacy of the safety procedures employed during the use of this product.

**End of Document**