

# SAFETY DATA SHEET

Version 6.6  
Revision Date 08/05/2024  
Print Date 08/06/2024

---

## SECTION 1: Identification of the substance/mixture and of the company/undertaking

### 1.1 Product identifiers

Product name : E-Cadherin Antibody (DECMA-1)  
Product Number : NB120-11512  
Brand : Novus Biologicals

### 1.2 Relevant identified uses of the substance or mixture and uses advised against

Identified uses : Laboratory chemicals, Synthesis of substances  
Uses advised against : The product is being supplied under the TSCA R&D Exemption (40 CFR Section 720.36). It is the recipient's responsibility to comply with the requirements of the R&D exemption. The product may not be used for a non-exempt commercial purpose under TSCA unless appropriate consent is granted in writing by Novus Biologicals.

### 1.3 Details of the supplier of the safety data sheet

Company : Novus Biologicals, LLC  
10771 E Easter Ave  
Centennial, CO 80112, USA  
Telephone : +1 888-506-6887  
Fax : +1 303-730-1966

### 1.4 Emergency telephone

Emergency Phone # : 888-506-6887

---

## SECTION 2: Hazards identification

### 2.1 Classification of the substance or mixture

Not a hazardous substance or mixture.

### 2.2 GHS Label elements, including precautionary statements

Not a hazardous substance or mixture.

### 2.3 Hazards not otherwise classified (HNOC) or not covered by GHS - none

---

**SECTION 3: Composition/information on ingredients****3.2 Mixtures**

Synonyms : Anti-E-Cadherin  
Anti-LCAM

| Component           | Classification        | Concentration   |
|---------------------|-----------------------|---|
| <b>sodium azide</b> |                       |   |
| CAS-No.             | 26628-22-8            | Acute Tox. 2; Acute Tox. 1; STOT RE 2; Aquatic Acute 1; Aquatic Chronic 1; H300, H330, H310, H373, H400, H410<br>M-Factor - Aquatic Acute: 1<br>M-Factor - Aquatic Chronic: 1 |
| EC-No.              | 247-852-1             |   |
| Index-No.           | 011-004-00-7          |   |
| Registration number | 01-2119457019-37-XXXX |   |
|                     |                       |   |
|                     |                       | >= 0.1 - < 1 %  |

For the full text of the H-Statements mentioned in this Section, see Section 16.

---

**SECTION 4: First aid measures****4.1 Description of first-aid measures**

No data available

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

The most important known symptoms and effects are described in the labelling (see section 2.2) and/or in section 11

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

No data available

---

**SECTION 5: Firefighting measures****5.1 Extinguishing media**

No data available

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

Nature of decomposition products not known.  
Not combustible.

**5.3 Advice for firefighters**

No data available

**5.4 Further information**

No data available

---

**SECTION 6: Accidental release measures****6.1 Personal precautions, protective equipment and emergency procedures**

For personal protection see section 8.

**6.2 Environmental precautions**

No data available

**6.3 Methods and materials for containment and cleaning up**

No data available

**6.4 Reference to other sections**

For disposal see section 13.

---

**SECTION 7: Handling and storage****7.1 Precautions for safe handling**

For precautions see section 2.2.

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

No data available

**7.3 Specific end use(s)**

Apart from the uses mentioned in section 1.2 no other specific uses are stipulated

---

**SECTION 8: Exposure controls/personal protection****8.1 Control parameters****Ingredients with workplace control parameters**

| Component    | CAS-No.    | Value                                  | Control parameters               | Basis   |
|--------------|------------|--|----------------------------------|---|
| sodium azide | 26628-22-8 | C                                      | 0.29 mg/m <sup>3</sup>           | USA. ACGIH Threshold Limit Values (TLV)   |
|              | Remarks    | Not classifiable as a human carcinogen |                                  |   |
|              |            | C                                      | 0.11 ppm                         | USA. ACGIH Threshold Limit Values (TLV)   |
|              |            | Not classifiable as a human carcinogen |                                  |   |
|              |            | C                                      | 0.1 ppm                          | USA. NIOSH Recommended Exposure Limits  |
|              |            | Potential for dermal absorption        |                                  |   |
|              |            | C                                      | 0.3 mg/m <sup>3</sup>            | USA. NIOSH Recommended Exposure Limits  |
|              |            | Potential for dermal absorption        |                                  |   |
|              |            | C                                      | 0.1 ppm<br>0.3 mg/m <sup>3</sup> | California permissible exposure limits for chemical contaminants (Title 8, Article 107) |
|              |            | Skin                                   |                                  |   |

## 8.2 Exposure controls

No data available

---

## SECTION 9: Physical and chemical properties

### 9.1 Information on basic physical and chemical properties

- |   |                              |
|---|------------------------------|
| a) Appearance                                   | Form: liquid                 |
| b) Odor   | No data available            |
| c) Odor Threshold                               | No data available            |
| d) pH   | No data available            |
| e) Melting point/freezing point                 | No data available            |
| f) Initial boiling point and boiling range      | No data available            |
| g) Flash point                                  | ( )No data available         |
| h) Evaporation rate                             | No data available            |
| i) Flammability (solid, gas)                    | No data available            |
| j) Upper/lower flammability or explosive limits | No data available            |
| k) Vapor pressure                               | No data available            |
| l) Vapor density                                | No data available            |
| m) Density                                      | No data available            |
| Relative density                                | No data available            |
| n) Water solubility                             | soluble                      |
| o) Partition coefficient: n-octanol/water       | No data available            |
| p) Autoignition temperature                     | Not applicable               |
| q) Decomposition temperature                    | No data available            |
| r) Viscosity                                    | No data available            |
| s) Explosive properties                         | Not classified as explosive. |
| t) Oxidizing properties                         | none                         |

### 9.2 Other safety information

No data available

---

## **SECTION 10: Stability and reactivity**

### **10.1 Reactivity**

No data available

### **10.2 Chemical stability**

No data available

### **10.3 Possibility of hazardous reactions**

No data available

### **10.4 Conditions to avoid**

No data available

### **10.5 Incompatible materials**

Strong oxidizing agents

### **10.6 Hazardous decomposition products**

In the event of fire: see section 5

---

## **SECTION 11: Toxicological information**

### **11.1 Information on toxicological effects**

#### **Mixture**

#### **Acute toxicity**

Acute toxicity estimate Oral - > 5,000 mg/kg

(Calculation method)

Acute toxicity estimate Inhalation - 4 h - 50.01 mg/l - dust/mist(Calculation method)

Acute toxicity estimate Dermal - > 5,000 mg/kg

(Calculation method)

#### **Skin corrosion/irritation**

No data available

#### **Serious eye damage/eye irritation**

No data available

#### **Respiratory or skin sensitization**

No data available

#### **Germ cell mutagenicity**

No data available

#### **Carcinogenicity**

IARC: No ingredient of this product present at levels greater than or equal to 0.1% is identified as probable, possible or confirmed human carcinogen by IARC.

NTP: No ingredient of this product present at levels greater than or equal to 0.1% is identified as a known or anticipated carcinogen by NTP.

OSHA: No component of this product present at levels greater than or equal to 0.1% is on OSHA's list of regulated carcinogens.

#### **Reproductive toxicity**

No data available

**Specific target organ toxicity - single exposure**

No data available

**Specific target organ toxicity - repeated exposure**

No data available

**Aspiration hazard**

No data available

**11.2 Additional Information**

To the best of our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated.

Hazardous properties cannot be excluded but are unlikely when the product is handled appropriately.

**Components**

**sodium azide**

**Acute toxicity**

LD50 Oral - Rat - 27 mg/kg

Remarks: (RTECS)

LC50 Inhalation - Rat - male and female - 4 h - 0.054 - 0.52 mg/l - dust/mist (US-EPA)

LD50 Dermal - Rabbit - 20 mg/kg

Remarks: (RTECS)

No data available

**Skin corrosion/irritation**

Skin - In vitro study

Result: No skin irritation

(OECD Test Guideline 439)

**Serious eye damage/eye irritation**

Eyes - Bovine cornea

Result: No eye irritation - 4 h

(OECD Test Guideline 437)

**Respiratory or skin sensitization**

Local lymph node assay (LLNA) - Mouse

Result: negative  
(OECD Test Guideline 429)

**Germ cell mutagenicity**

Test Type: Mutagenicity (mammal cell test): chromosome aberration.

Test system: Chinese hamster ovary cells

Result: negative

Test Type: unscheduled DNA synthesis assay

Test system: Chinese hamster lung cells

Result: negative

Test Type: sister chromatid exchange assay

Test system: Chinese hamster ovary cells

Result: negative

**Carcinogenicity**

No data available

**Reproductive toxicity**

No data available

**Specific target organ toxicity - single exposure**

No data available

**Specific target organ toxicity - repeated exposure**

Oral - May cause damage to organs through prolonged or repeated exposure.

- Brain

**Aspiration hazard**

No data available

---

**SECTION 12: Ecological information**

**12.1 Toxicity**

**Mixture**

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**

PBT/vPvB assessment not available as chemical safety assessment not required/not conducted

**12.6 Endocrine disrupting properties**

No data available

**12.7 Other adverse effects**

No data available

## Components

### sodium azide

|                      |   |
|----------------------|---|
| Toxicity to fish     | flow-through test LC50 - Oncorhynchus mykiss (rainbow trout)<br>- 2.75 mg/l - 96 h<br>(OECD Test Guideline 203) |
| Toxicity to algae    | static test ErC50 - Pseudokirchneriella subcapitata - 0.35 mg/l<br>- 96 h<br>(OECD Test Guideline 201)          |
| Toxicity to bacteria |   |

---

## SECTION 13: Disposal considerations

### 13.1 Waste treatment methods

No data available

---

## SECTION 14: Transport information

### DOT (US)

Not dangerous goods

### IMDG

Not dangerous goods

### IATA

Not dangerous goods

### Further information

Not classified as dangerous in the meaning of transport regulations.

---

## SECTION 15: Regulatory information

### CERCLA Reportable Quantity

| Components   | CAS-No.    | Component RQ (lbs) | Calculated product RQ (lbs) |
|--------------|------------|--------------------|-----------------------------|
| sodium azide | 26628-22-8 | 1000               |                             |

### SARA 304 Extremely Hazardous Substances Reportable Quantity

| Components   | CAS-No.    | Component RQ (lbs) | Calculated product RQ (lbs) |
|--------------|------------|--------------------|-----------------------------|
| sodium azide | 26628-22-8 | 1000               |                             |



## **SARA 302 Extremely Hazardous Substances Threshold Planning Quantity**

This material does not contain any components with a section 302 EHS TPQ.

**SARA 311/312 Hazards** : No SARA Hazards

**SARA 313** : This material does not contain any chemical components with known CAS numbers that exceed the threshold (De Minimis) reporting levels established by SARA Title III, Section 313.

### **US State Regulations**

#### **Massachusetts Right To Know**

sodium azide 26628-22-8

#### **Massachusetts Right To Know**

water 7732-18-5  
sodium azide 26628-22-8

#### **Pennsylvania Right To Know**

water 7732-18-5  
disodium hydrogen orthophosphate 7558-79-4  
sodium azide 26628-22-8

#### **Pennsylvania Right To Know**

sodium azide 26628-22-8

#### **Maine Chemicals of High Concern**

water 7732-18-5  
sodium chloride 7647-14-5  
disodium hydrogen orthophosphate 7558-79-4  
sodium dihydrogen phosphate 7558-80-7

#### **Vermont Chemicals of High Concern**

water 7732-18-5  
sodium chloride 7647-14-5  
disodium hydrogen orthophosphate 7558-79-4  
sodium dihydrogen phosphate 7558-80-7

#### **Washington Chemicals of High Concern**

water 7732-18-5  
sodium chloride 7647-14-5  
disodium hydrogen orthophosphate 7558-79-4  
sodium dihydrogen phosphate 7558-80-7

#### **New Jersey Right To Know**

water 7732-18-5

### **The ingredients of this product are reported in the following inventories:**

TSCA : Product contains substance(s) not listed on TSCA inventory.

**TSCA list**

No substances are subject to a Significant New Use Rule.

No substances are subject to TSCA 12(b) export notification requirements.

---

**SECTION 16: Other information**

The branding on the header and/or footer of this document may temporarily not visually match the product purchased as we transition our branding. However, all of the information in the document regarding the product remains unchanged and matches the product ordered.

Version: 6.6

Revision Date: 08/05/2024

Print Date: 08/06/2024