

Date of Issue: 05/29/20

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Section 1. Identification of the Substance/Mixture and of the Company/Undertaking

1.1 Product Identifier

Substance Name/Product Name: QuantTag Biotin Quantitation Kit – Reagent 3

Product Identifier/Catalog No.: BDK-2000

CAS Number: Not Applicable/Chemical Mixture

EC Number: Not Applicable/Chemical Mixture

REACH Registration Number: Not Applicable/Chemical Mixture

1.2 Relevant identified uses and uses advised against

Relevant identified uses: Research Grade Use Only, For professional Users Only

Uses advised against: None

Reason why uses advised against: Not Applicable

1.3 Details of Safety Data Sheet Supplier

Manufacturer/Supplier USA:

Vector Laboratories, Inc.

30 Ingold Road

Burlingame, CA 94010

United States

Tel: (650) 697-3600 Fax: (650) 697-0339

Email: vector@vectorlabs.com

Supplier United Kingdom:

Vector Laboratories Ltd.

3 Accent Park

Bakewell Road, Orton Southgate

Peterborough, PE2 6XS, United Kingdom

Tel: (01733) 237999 Fax: (01733) 237119

Email: vector@vectorlabs.co.uk

1.4 Emergency Telephone Number

Emergency Telephone Number: (650) 697-3600

Opening hours: 5 Days a Week – 9 am to 4 pm PST/PDT

Section 2. Hazards Identification

GHS Classification:

Flammable Liquids - Category 2, H225

Acute Toxicity, Oral - Category 4, H302

Acute Toxicity, Inhalation - Category 4, H332

Acute Toxicity, Dermal - Category 4, H312

Eye irritation - Category 2A, H319

GHS Label:



Signal Word:

Danger

Hazard Statement(s):

H225 Highly flammable liquid and vapor.

H302 + H312 + H332 Harmful if swallowed, in contact with skin or if inhaled.

H319 Causes serious eye irritation

Date of Issue: 05/29/20

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Precautionary Statement(s):

Keep away from heat/sparks/open flames/hot surfaces. - NO SMOKING. Keep container tightly closed. Use explosion-proof electrical/ventilating/lighting/equipment. Wear protective gloves/protective clothing/eye protection/face protection. Do not breathe dust/fume/gas/mist/vapors/spray. Wash skin thoroughly after handling. IF SWALLOWED: Call a POISON CENTER or doctor/physician if you feel unwell. Rinse mouth. IF ON SKIN: Take off immediately all contaminated clothing. Rinse skin with water/shower. IF INHALED: Remove person to fresh air and keep comfortable for breathing. Call a POISON CENTER or doctor/physician if you feel unwell. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. If eye irritation persists: Get medical advice/attention.

Section 3. Composition/Information on Ingredients

Component	Identifier	Classification	% (Weight)
Acetonitrile	CAS # 75-05-8 EC # 200-835-2	Flammable Liquids - Category 2, H225 Acute Toxicity, Oral - Category 4, H302 Acute Toxicity, Inhalation - Category 4, H332 Acute Toxicity, Dermal - Category 4, H312 Eye irritation - Category 2A, H319	>90%

Section 4. First Aid Measures

Inhalation: Remove to fresh air. Obtain medical attention.

Skin contact: Wash thoroughly with water for at least 15 minutes. Obtain medical attention.

Eye contact: Wash thoroughly with water for at least 15 minutes. Obtain medical attention.

Ingestion: Do not induce vomiting. If person is conscious, wash out mouth with water. Obtain medical attention.

Section 5. Fire Fighting Measures

Suitable extinguishing media: Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.

Special protective equipment: Use respirator. Toxic gases can be evolved.

Section 6. Accidental Release Measures

Wear protective clothing to prevent contact with eyes and skin. Ensure adequate ventilation. Collect on absorbent material and dispose of according to federal, state and local environmental regulations.

Section 7. Handling and Storage

Handling: Use only with adequate ventilation. Wear eye protection and compatible chemical-resistant gloves.

Storage: Keep container tightly closed and store in a refrigerator.

Date of Issue: 05/29/20

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Section 8. Exposure Controls/Personal Protection**Exposure limits:** Use enclosures**Engineering controls:**

Safety shower and eye bath. Mechanical exhaust required.

Hand protection - Compatible chemical-resistant gloves.**Eye protection** - Chemical safety goggles.**Skin protection** - Lab coat.**Section 9. Physical and Chemical Properties****Appearance:** Colorless liquid**Flash point:** 5.6 °C**Odor:** None**Autoflammability:** Not determined**pH:** Not determined**Vapor pressure:** Not determined**Boiling point:** 81.6 °C**Relative density:** Not determined**Melting point:** -44 °C**Solubility:** Not determined**Section 10. Stability and Reactivity****Chemical stability:** Stable under recommended storage conditions.**Conditions to avoid:** Heat, sparks, open flame.**Materials to avoid:** Strong oxidizing agents.**Hazardous reactions/decomposition products:** Carbon monoxide, carbon dioxide, nitrogen oxides, hydrogen.**Section 11. Toxicological Information****Information on toxicological effects:****Acute effects:** No data available.**Serious Eye Damage/Eye Irritation:** Irritating to eyes.**Reproductive Toxicity:** No data available.**STOT - SE:** The substance or mixture is not classified as specific target organ toxicant.**Acute effects:** Harmful if inhaled. May cause respiratory tract irritation. Harmful if absorbed through skin. May cause skin irritation. Causes eye irritation. Harmful if swallowed.**Chronic effects:** Not determined.

Date of Issue: 05/29/2

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

RTECS#: AL7700000

Toxicity data: orl-rat LD50: 3800 mg/kg
skn-rbt LD50: 1250 mg/kg
orl-gpg LD 50: 177 mg/kg

Only selected RTECS data is shown.

Section 12. Ecological Information

Avoid release to the environment.

Section 13. Disposal Considerations

Observe all federal, state and local environmental disposal regulations.

Section 14. Transport Information

DOT (US): Excepted Quantities: Packing group: II
UN number: 1648 Class: 3
Proper shipping name Acetonitrile
Reportable Quantity (RQ): 5000 lbs.

DOT (US): Excepted Quantities: Packing group: II
UN number: 1648 Class: 3
Proper shipping name Acetonitrile

IMDG: Excepted Quantities: Packing group: II
UN number: 1648 Class: 3
Proper shipping name Acetonitrile

IATA: Excepted Quantities: Packing group: II
UN number: 1648 Class: 3
Proper shipping name Acetonitrile

Section 15. Regulatory Information

SARA:

SARA 302 - No chemicals in this material are subject to the reporting requirements.

SARA 313 - The following components are subject to reporting levels: Acetonitrile, CAS-No. 75-05-8

SARA 311/312 Hazards - Fire Hazard, Acute Health Hazard

Risk phrases:

R11 - Highly flammable.

R20/21/22 - Harmful by inhalation, in contact with skin and if swallowed

R36 - Irritating to eyes

Date of Issue: 05/29/2

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Safety phrases:

S16 - Keep away from sources of ignition - No Smoking

S36/37 - Wear suitable protective clothing and gloves.

California Prop. 65 components:

This product does not contain any chemical known to the State of California to cause cancer, birth defects, or other reproductive harm

Section 16. Other Information

The information contained in this Material Safety Data is believed to be accurate but it is the responsibility of the user to determine the applicability of these data to the formulation of necessary safety precautions. Vector Laboratories, Inc. shall not be held responsible for any damage resulting from the use of the above product or the information contained in this Material Safety Data Sheet.

Date of Issue: 05/29/2

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Section 1. Product and Company Identification

Product Name: QuantTag Biotin Quantitation Kit – Other Kit components
(Reagent 1, Reagent 2, & Biotin Standard Solution)

Catalog No.: BDK-2000

Manufacturer/Supplier USA:

Vector Laboratories, Inc.
30 Ingold Road
Burlingame, CA 94010
United States
Tel: (650) 697-3600 Fax: (650) 697-0339
Email: vector@vectorlabs.com

Supplier United Kingdom:

Vector Laboratories Ltd.
3 Accent Park
Bakewell Road, Orton Southgate
Peterborough, PE2 6XS, United Kingdom
Tel: (01733) 237999 Fax: (01733) 237119
Email: vector@vectorlabs.co.uk

Section 2. Hazards Identification

GHS Classification:

These components have been classified as non-hazardous based on the physical and/or chemical nature and/or concentration of ingredients.

GHS Label: Non-hazardous substance or mixture.

Hazards not otherwise classified (HNOC) or not covered by GHS: None

Section 3. Composition/Information on Ingredients

No ingredients are hazardous according to OSHA criteria.
No components need to be disclosed according to the applicable regulations.

Section 4. First Aid Measures

Inhalation: Remove to fresh air.
Skin contact: Wash thoroughly with soap and water.
Eye Contact: Flush eyes with water as a precaution.
Ingestion: Do not induce vomiting. Wash out mouth with water.

Section 5. Fire Fighting Measures

Suitable extinguishing media: Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.
Special protective equipment: Use respirator for fire fighting if necessary.

Section 6. Accidental Release Measures

Date of Issue: 05/29/2

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Wear protective clothing to prevent contact with eyes and skin. Ensure adequate ventilation. Collect on absorbent material and dispose of according to federal, state and local environmental regulations.

Section 7. Handling and Storage

Handling: Use only with adequate ventilation. Wear eye protection and compatible chemical-resistant gloves. Handle in accordance with good industrial hygiene and safety practice.

Storage: Keep container tightly closed and store according to product label.

Section 8. Exposure Controls/Personal Protection

Exposure limits: Contains no substances with occupational exposure limit values.

Engineering controls: Use enclosures, local exhaust ventilation or other engineering controls to keep airborne levels below recommended exposure limits.

Personal protection:

Hand protection - Compatible chemical-resistant gloves.

Eye protection - Chemical safety goggles.

Skin protection - Lab coat.

Section 9. Physical and Chemical Properties

Appearance: Liquid

Odor: None

pH: Not determined

Boiling point: Not determined.

Melting point: Not determined.

Flash Point: Not determined.

Autoflammability: Not determined.

Vapor pressure: Not determined.

Relative density: Not determined.

Solubility: Not applicable.

Section 10. Stability and Reactivity

Chemical stability: Stable under recommended storage conditions.

Conditions to avoid: pH extremes.

Materials to avoid: Not applicable.

Hazardous reactions/decomposition products: Not determined.

Section 11. Toxicological Information

Acute toxicity: No data available.

Chronic toxicity: No data available.

Section 12. Ecological Information

No data available.

Section 13. Disposal Considerations

Observe all federal, state and local environmental regulations.

Date of Issue: 05/29/2

Version 1.0

Previous Date of Issue: 04/15/15

Previous Version: N/A

Section 14. Transport Information**DOT (US):** Not dangerous goods.**RID/ADR:** Not regulated.**IMDG:** Not dangerous goods.**IATA:** Not dangerous goods.**Section 15. Regulatory Information****Substances of Very High Concern:** None**SARA:**

SARA 302 - No chemicals in this material are subject to the reporting requirements.

SARA 313 - No chemicals in this material with known CAS #s are subject to the reporting requirements.

SARA 311/312 Hazards - No SARA Hazards.

California Prop. 65 components:

This product does not contain any chemicals known to the state of California to cause cancer, birth defects or other reproductive harm.

Section 16. Other Information

The information contained in this Safety Data Sheet is believed to be accurate, but it is the responsibility of the user to determine the applicability of this data to necessary safety precautions. Vector Laboratories, Inc. shall not be held responsible for any damage resulting from the use of the above product or the information contained in this Safety Data Sheet.

For Research Grade Use Only.