

SAFETY DATA SHEET

Date of Issue: August 23, 2018

Section 1. Product and Company Identification

Product Name: VECTASHIELD® Vibrance™ Antifade Mounting Medium

Catalog No.: H-1700

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

For Research Grade Use Only.

Section 2. Hazards Identification

GHS Classification:

This product has 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.



SAFETY DATA SHEET

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.

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. Flash Point: Not determined.

Odor: None.

pH: 8.1

Vapor pressure: Not determined.

Boiling point: Not determined.

Relative density: Not determined.

Melting point: 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.

H-1700



SAFETY DATA SHEET

Section 13. Disposal Considerations

Observe all federal, state and local environmental regulations.

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

Date of Issue or Revision: August 23, 2018

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.

H-1700