



MATERIAL SAFETY DATA SHEET

Date of Issue: May 6, 2008

1. Identification of the substance/preparation and of the company/undertaking

Product Name: VECTASTAIN® ABC-AmP™ KITS
Catalog No.: AK-6000, 6401, 6402, 6601, 6602

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

EU 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

2. Hazards identification

Highly Toxic (USA) Very Toxic (EU). Dangerous for the environment. Very toxic by inhalation, in contact with skin and if swallowed. Readily absorbed through skin. Contact with acids liberates very toxic gas. Very toxic to aquatic organism organisms, may cause long-term adverse effects in the aquatic environment.

3. Composition/information on ingredients

Chemical characterization: Mixture of substance(s) listed below with non-hazardous additions

Dangerous chemical	CAS#	EC#	% by weight
sodium azide	26628-22-8	247-852-1	0.08%

Classification

T+-N; R28-32-50/53

4. First-aid measures

Inhalation: Remove to fresh air. Obtain medical attention.

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

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

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

5. Fire-fighting measures

Suitable extinguishing media: Water, foam, carbon dioxide

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

6. Accidental release measures

Wear protective clothing to prevent contact with eyes and skin. Ensure adequate ventilation. Flush to sewer, if allowed or collect on absorbent material and dispose of according to local and regional regulations.



7. Handling and storage

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

Storage: Store in a refrigerator.

8. Exposure controls/personal protection

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

Exposure limit: TLV 0.29 mg/m³ from ACGIH. Consult local authorities for acceptable exposure limits.

Personal protection:

Hand protection Compatible chemical-resistant gloves

Eye protection Chemical safety goggles

Skin protection Lab coat

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

10. Stability and reactivity

Conditions to avoid: acidic pH

Materials to avoid: strong oxidizers

Hazardous decomposition products: carbon monoxide, carbon dioxide, nitrogen oxides.

11. Toxicological information

Acute effects: May cause skin irritation, eye irritation, mucus membrane irritation. May be fatal if swallowed, inhaled or absorbed through the skin. Exposure can cause nausea, headache and vomiting.

Chronic effects: May alter genetic material

RTECS#: VY8050000

Toxicity data: orl-hmn TDLo: 710 µg/kg

orl-rat LD50: 27 mg/kg

ivn-mus LD50: 19 mg/kg

skn-rbt LD50: 20 mg/kg

Only selected RTECS data is shown. See actual entry for complete information.

(continued on next page)



12. Ecological information

Data not yet available.

13. Disposal considerations

Observe all local or regional regulations.

14. Transport information

Data not yet available.

15. Regulatory information

Hazard symbols:

T+ Very toxic

N Dangerous for the environment

Risk phrases:

R28 Very toxic if swallowed

R32 Contact with acids liberates very toxic gas

R50/53 Very toxic to aquatic organism organisms, may cause long-term adverse effects in the aquatic environment

Safety phrases:

S28 After contact with skin, wash immediately with plenty of soap-suds

S45 In case of accident or if you feel unwell, seek medical advice immediately (show the label where possible)

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.