

TREVIGEN®

MATERIAL SAFETY DATA SHEET

CHEMICAL IDENTIFICATION

PRODUCT NAME: Cultrex® Bovine Collagen I (in 20 mM Acetic Acid)

CATALOG NUMBERS: 3442-005-01; 3442-050-01

COMPOSITION/INFORMATION ON INGREDIENTS

Bovine Collagen I	CAS#: N/A	SARA 313: No
Acetic Acid (< 1%)	CAS#: 64-19-7	SARA 313: No

HAZARDS IDENTIFICATION

Emergency Overview:

OSHA Hazards: Target Organ Effect, Harmful by skin absorption, Skin sensitizer, Corrosive

Target Organs: Teeth, Kidney

GHS Classification:

Skin irritation (Category 2)
Eye irritation (Category 2A)
Skin sensitization (Category 1)

Hazard statements:

H315 Causes skin irritation
H317 May cause an allergic skin reaction
H319 Causes serious eye irritation

Precautionary Statements:

P280 Wear protective gloves
P305 + P351 + P338 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

HMIS Classification:	Health: 3	Flammability: 0	Reactivity: 0
NFPA Rating:	Health: 3	Flammability: 0	Reactivity: 0

FIRST AID MEASURES

Eye exposure: In case of contact, flush eyes with plenty of water for 15 minutes. Continue rinsing eyes during transport to hospital. Consult a physician.

Oral exposure: Never give anything by mouth to an unconscious person. Rinse mouth with water. Consult a physician.

Dermal exposure: In case of contact, wash with plenty of soap and water. Consult a physician.

Inhalation exposure: If inhaled, remove to fresh air. If breathing with difficulty, give oxygen. If not breathing, give artificial respiration.

FIRE FIGHTING MEASURES

FLAMMABILITY: Not flammable or combustible.

EXTINGUISHING MEDIA

Suitable: Water spray, Carbon dioxide, dry chemical powder, or alcohol-resistant foam.

Hazardous combustion products: Carbon oxides

PROCEDURE(S) OF PERSONAL PRECAUTION(S)

Avoid contact with eyes and skin. Avoid inhalation of vapor or mist. Wear laboratory coat, safety goggles, rubber boots, and laboratory rubber gloves. Wash contaminated clothing thoroughly after handling and before reuse.

METHODS FOR CLEANING UP

Absorb on sand or vermiculite and place in closed containers for disposal as hazardous waste. Ventilate area and wash spill site after material pickup is complete.

CONTINUED ON NEXT PAGE

STORAGE

Store at 4 °C. Opened containers must be carefully resealed and kept upright to prevent leakage.

STABILITY

Chemically stable under recommended storage conditions.

Materials to Avoid: Oxidizing agents, soluble carbonates and phosphates, hydroxides, metals, peroxides, permanganates, e.g. potassium permanganate, amines, alcohols.

ROUTE OF EXPOSURE

Inhalation: May be harmful; material is destructive to the tissue of the mucous membranes and upper respiratory tract.
 Skin: Causes skin burns.
 Eyes: Causes eye burns.
 Ingestion: May be harmful if swallowed.

When used according to the instructions for use, this product does not have harmful effects according to our experience and to our knowledge.

DISPOSAL

Observe all federal, state, and local environmental regulations. Offer surplus and non-recyclable solutions to a licensed disposal company. Contact a licensed professional waste disposal service to dispose of this material. Dissolve or mix the material with a combustible solvent and burn in a chemical incinerator equipped with an afterburner and scrubber.

SHIPPING

DOT: Not dangerous goods
 IMDG: Not dangerous goods
 IATA: Not dangerous goods

REGULATORY INFORMATION

OSHA Hazards: Target Organ Effect, Harmful by skin absorption, skin sensitizer, corrosive

SARA 302: No chemicals in this material are subject to the reporting requirements of SARA Title III, Section 302

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.

SARA 311/312: Acute Health Hazard, Chronic Health Hazard

Massachusetts Right to Know Components

Acetic Acid

CAS-No.

64-19-7

Revision Date

1993-04-24

Pennsylvania Right to Know Components

Acetic Acid

CAS-No.

64-19-7

Revision Date

1993-04-24

New Jersey Right to Know Components

Acetic Acid

CAS-No.

64-19-7

Revision Date

1993-04-24

California Prop. 65 Components

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

CONTINUED ON NEXT PAGE

OTHER INFORMATION

The above information is believed to be correct, however, the physical, chemical, and toxicological properties of the product may not yet have been fully characterized. Trevigen recommends the use of gloves, lab coats and eye protection while using or handling this product. Trevigen assumes no liability for damage resulting from handling or contact with the above product. This MSDS does not represent any guarantee of the properties of the product.

Trevigen, Inc.

8405 Helgerman Ct.

Gaithersburg, MD 20877

Tel: 1-800-873-8443 • 301-216-2800

Fax: 301-560-4973

e-mail: info@trevigen.comwww.trevigen.com