

## MATERIAL SAFETY DATA SHEET

Date: March 21, 2016

## 1. PRODUCT AND COMPANY IDENTIFICATION

**Product Name:** Annexin V, FITC conjugate (FITC Annexin V)  
**Catalog Number:** 29001  
**Unit Size:** 0.5mL  
**Manufacturer/Supplier:** Biotium, Inc.  
 46117 Landing Parkway, Fremont, CA 94538, USA  
 Phone: 1-510-265-1027, Fax: 1-510-265-1352  
 Web: <http://www.biotium.com>

## 1. HAZARDS IDENTIFICATION

**Classification according to Regulation (EC) No 1272/2008[CLP]**

None

**Classification according to Directive 1999/45/EC**

None

**Labelling according to Regulation (EC) No 1272/2008[CLP]****Hazard pictogram**

None

**Signal word** None**Hazard statements** None**Precautionary statements** None**HMIS Classification**

Health hazard: 0

Flammability: 0

Physical hazards: 0

**NFPA Rating**

Health hazard: 0

Fire: 0

Reactivity Hazard: 0

## 2. COMPOSITION/INFORMATION ON INGREDIENTS

## Liquid Format CF dye Annexin V conjugates

Name	CAS No.	EC No.	Index No.	Weight %	Classification according to 67/548/EEC	Classification according to regulation (EC)No1278/2008
FITC Annexin V	-	-	-	<0.05%	-	-
Bovine Serum Albumin	9048-46-8	232-936-2	-	<0.5%	-	-
Tris Base	77-86-1	201-064-4	-	<0.2%	-	-
EDTA	60-00-4	200-449-4	607-429-00-8	<0.1%	Xi; R36	-Eye Irrit. 2

						H319
Sodium Azide	26628-22-8	247-852-1	011-004-00-7	<0.1%	T+= very toxic, R28,R32,R50-53	Acute Tox.2, H300,H400,H410

---

### 3. FIRST-AID MEASURES

**General advice**

Consult a physician. Show this safety data sheet to the doctor in attendance. Move out of dangerous area.

**If inhaled**

If breathed in, move person into fresh air. If not breathing, give artificial respiration. Consult a physician.

**In case of skin contact**

Wash off with soap and plenty of water. Consult a physician.

**In case of eye contact**

Flush eyes with water as a precaution.

**If swallowed**

Do NOT induce vomiting. Never give anything by mouth to an unconscious person. Rinse mouth with water. Consult a physician.

---

### 4. FIREFIGHTING MEASURES

**Suitable extinguishing media**

carbon dioxide, dry chemical extinguishers, foam extinguishers or water.

**Special protective equipment for firefighters**

Wear self contained breathing apparatus for firefighting if necessary.

---

### 5. ACCIDENTAL RELEASE MEASURES

**Personal precautions**

Avoid breathing vapors, mist or gas. Remove all sources of ignition.

**Environmental precautions**

Prevent further leakage or spillage if safe to do so. Do not let product enter drains.

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

Contain spillage. Soak up spilled substance with inert absorbent material. . Keep in suitable, closed containers for disposal.

---

### 6. HANDLING AND STORAGE

**Precautions for safe handling**

Avoid inhalation of vapor or mist.

Avoid direct contact with substance.

**Conditions for safe storage**

Keep container tightly closed in a dry and well-ventilated place.

Store at 4 °C.

## 7. EXPOSURE CONTROLS/PERSONAL PROTECTION

### Components with workplace control parameters

Substance Sodium Azide

CAS no. 26628-22-8

country	Austria	Belgium	Denmark	European Union	France	Germany
Limit value, 8hours	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.2mg/m <sup>3</sup>
Limit value, short term	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	0.2mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	0.4mg/m <sup>3</sup>

country	Hungary	Italy	Poland	Spain	Sweden	Netherlands	United Kingdom
Limit value, 8hours	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>	-	0.1mg/m <sup>3</sup>	0.1mg/m <sup>3</sup>
Limit value, short term	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	-	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>

country	USA-NIOSH	USA-OSHA	Australia	Canada	Japan	South Korea
Limit value, 8hours	-	-	-	-	-	-
Limit value, short term	0.3mg/m <sup>3</sup>	-	0.3mg/m <sup>3</sup>	0.3mg/m <sup>3</sup>	-	-

### Personal protective equipment

#### Hand protection

Handle with gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.

#### Respiratory protection

Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

#### Eye protection

Safety glasses with side-shields conforming to EN166 Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU).

#### Hygiene measures

Handle in accordance with good industrial hygiene and safety practice. Wash hands before breaks and at the end of workday.

## 8. PHYSICAL AND CHEMICAL PROPERTIES

Chemical Name	FITC Annexin V
Appearance	Liquid
Odor	No data available
Odor threshold	No data available
pH	7-8
Melting point/freezing point	No data available
Boiling point	No data available
Flash point	No data available
Evaporate rate	No data available
Flammability	No data available

Explosive limits	No data available
Vapor pressure	No data available
Vapor density	No data available
Relative density	No data available
Solubility	Water soluble
Partition coefficient:n-octanol/water	No data available
Auto-ignition temperature	No data available
Decomposition temperature	No data available
Viscosity	No data available
Explosive properties	No data available
Oxidising properties	No data available

---

## 9. STABILITY AND REACTIVITY

**Chemical stability**

Stable under recommended storage conditions.

**Possibility of hazardous reactions**

No data available

**Conditions to avoid**

Heat, flames and sparks.

**Materials to avoid**

No data available

**Hazardous decomposition products**

Hazardous decomposition products formed under fire conditions. - Carbon oxides

Other decomposition products - no data available

---

## 10. TOXICOLOGICAL INFORMATION

**Acute toxicity**

Chemical Name	LD50 (oral) rat	LD50 (dermal) rabbit	LC50 (inhalation)
Sodium Azide	27mg/kg	20mg/kg	No data available

**Other information on acute toxicity** no data available

**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 component of this product present at levels greater than or equal to 0.1% is identified as probable, possible or confirmed human carcinogen by IARC.

ACGIH: No component of this product present at levels greater than or equal to 0.1% is identified as a carcinogen or potential carcinogen by ACGIH.

NTP: No component 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 identified as a carcinogen or potential carcinogen by OSHA.

**Reproductive toxicity** no data available

**Specific target organ toxicity - single exposure (Globally Harmonized System)**

no data available

**Specific target organ toxicity - repeated exposure (Globally Harmonized System)**

no data available

**Aspiration hazard**

no data available

**Potential health effects**

Inhalation May be harmful if inhaled. May cause respiratory tract irritation.

Ingestion May be harmful if swallowed.

Skin May be harmful if absorbed through skin. May cause skin irritation.

Eyes May cause eye irritation.

**Additional Information**

RTECS: None

---

**11. ECOLOGICAL INFORMATION****Toxicity** no information available**Persistence and degradability** no information available**Biodegradation** no information available**Mobility in soil** no information available**Results of PBT and vPvB assessment** no information available**Other adverse effects** no information available**Additional information** no information available

---

**12. DISPOSAL CONSIDERATIONS**

Do not dispose product directly into sewage. Consult local state or national regulation for proper disposal.

---

**13. TRANSPORT INFORMATION****IATA, IMDG, DOT(US)** not dangerous good during transportation

UN number none

UN proper shipping name none

Transport hazard class none

Packing group none

Environmental hazards none

Special precaution for user none

---

**14. REGULATION INFORMATION****US Federal Regulations**

Us Toxic Substances Control Act(TSCA): Not listed

SARA 302: No chemicals were found .

SARA 313: No chemicals were found.

SARA 311/312 Hazards : No chemicals were found.

---

**15. OTHER INFORMATION**

Classification according to Regulation (EC) Nr. 1272/2008

Prepared by: Regulatory Department  
Biotium Inc.  
Version no. 4  
Reason for revision Application of CLP labeling and corresponding requirements

The information provided above is believed to be correct to our best knowledge, but does not purport to be all inclusive, and shall be used only as a guide. This material is sold for research purposes only and is not required to appear on the TSCA inventory. It is not intended for food, drug, household, agricultural or cosmetic use. Its use must be supervised by a technically qualified individual experienced in handling potentially hazardous chemicals. Biotium shall not be held liable for any damage resulting from handling or contact with the above product.