

SAFETY DATA SHEET

Section 1 – Identification

Product Name	FITC Annexin V Apoptosis Detection Kit with PI
Catalog No.	640914
Recommended Use	Research use only.
Company	BioLegend
Street Address	9727 Pacific Heights Blvd
City, State, Zip, Country	San Diego, CA 92121
Phone	+1-858-455-9588
Emergency Number	In case of a chemical emergency, spill, fire, or exposure, +1-858-455-9588 (7:00AM – 5:00PM PDT, M-F)

Section 2 – Hazards Identification

2.1 Hazard Classification

Not classified

2.2 GHS Label elements, including precautionary statements

Pictogram None

Signal Word None

Hazard Statement None

Precautionary Statement None

Section 3 – Composition/Information on Ingredients

Component	CAS	EINECS	EU Index	Concentration
Propidium Iodide	25535-16-4	N/A	N/A	<0.1%
Sodium Azide	26628-22-8	247-852-1	011-004-00-7	0.09%

Section 4 – First Aid Measures

4.1 Description of first aid measures

General information: Symptoms of poisoning may even occur after several hours; therefore provide medical observation for at least 48 hours after the accident.

After inhalation: Move to fresh air. If breathing is difficult/has stopped, give artificial respiration using respiratory protection. Do not use mouth-to-mouth method of resuscitation. Seek medical attention.

After skin contact: Immediately remove contaminated clothes, and wash with soap and copious amounts of water. Seek immediate medical attention.

After eye contact: Immediately flush eyes with plenty of water for at least 15 minutes, occasionally lifting the eye lids to ensure thorough rinsing. Seek medical attention.

After swallowing: Wash mouth out with water if person is conscious. DO NOT induce vomiting. Drink copious amounts of water, milk, or milk magnesia if conscious. Seek medical attention.

Symptoms: Contact may cause skin irritation. Contact may cause eye irritation.

Section 5 – Fire-Fighting Measures

5.1 Suitable extinguishing agents: CO₂, extinguishing powder or water spray. Fight larger fires with water spray or alcohol resistant foam.

5.2 Special hazards caused by the material, its products of combustion or resulting gases: During heating or in case of fire, poisonous gases may be produced.

5.3 Special protective equipment and precautions for fire-fighters: Wear protective clothing and self-contained breathing apparatus.

Hazardous combustion materials: No data available.

Section 6 – Accidental Release Measures

6.1 Personal precautions, protective equipment, and emergency procedures: Not required

6.2 Environment precautions: Prevent entry into waterways, drains, soil, and sewers. Use suitable absorbent material.

6.3 Measures for cleaning/collecting: Absorb with liquid-binding material (sand, diatomite, acid binders, universal binders). Do not allow to enter any water system.

6.4 Additional information:

See Section 7 for information on safe handling.

See Section 8 for information on personal protection equipment.

See Section 13 for disposal information.

Section 7 – Handling and Storage

7.1 Information for safe handling: No special measures required.

7.2 Conditions for safe storage, including any incompatibilities: Do not store together with oxidizing and acidic materials as well as heavy-metal compounds.

Section 8 – Exposure controls/personal protection

8.1 Exposure Limits

Sodium Azide

Cal/OSHA

PEL

Ceiling - 0.1 ppm, 0.3 mg/m³

ACGIH
TLV Ceiling – 0.29 mg/m³

NIOSH
REL Ceiling – 0.3 mg/m³

8.2 Exposure Controls

Engineering Controls

Use only with adequate (local exhaust) ventilation or fume hood.

Personal protective equipment

General protective and hygienic measures

Keep away from foodstuffs, beverages, and feed.

Wash hands, face, and exposed forearms/areas after handling.

Wash contaminated clothing before reusing.

Ensure eyewash stations and safety showers are in close proximity to workstation.

Breathing equipment: No required.

Protection of hands: Chemical resistant gloves.

Eye protection: Safety goggles.

Body protection: Protective work clothing (lab coat).

Section 9 – Physical and Chemical Properties

Appearance	Liquid, Colored, clear
Odor	Odorless
Odor Threshold	Not Available
pH	7.0-7.2
Melting point/freezing point	No Data Available
Boiling point	No Data Available
Flash point	No Data Available
Evaporation rate	No Data Available
Flammability	No Data Available
Upper explosion limit	No Data Available
Lower explosion limit	No Data Available
Vapor pressure	No Data Available
Vapor density	No Data Available
Relative density	No Data Available
Solubility	Soluble
Partition coefficient	No Data Available
Auto-ignition temperature	No Data Available
Decomposition temperature	No Data Available
Viscosity	No Data Available
Explosive Properties	No Data Available
Oxidizing Properties	No Data Available

Section 10 – Stability and Reactivity

10.1 Reactivity

No data available

10.2 Chemical stability

Stable under normal ambient and anticipated storage and handling conditions of temperature and pressure.

10.3 Possibility of hazardous reactions

No data available

10.4 Conditions to avoid

No data available

10.5 Incompatible materials

Acidic materials and metals.

10.6 Hazardous decomposition products

Nitrogen Oxides (NO_x)

10.5 Additional information

Sodium azide is present in this product. Contact with acidic solutions and metal compounds over time may form potentially explosive metal azides. Should any of this material be introduced into a sanitary sewer system, flush with copious amounts of water.

Section 11 – Toxicological Information

11.1 Information on toxicological effects

Routes of Entry	Ingestion, inhalation, skin and eye contact.
Acute Toxicity	Oral LD50 (Sodium Azide) 27 mg/kg (rat)
Skin Corrosion/Irritation	Contact may cause skin irritation
Serious eye damage/irritation	Contact may cause eye irritation
Respiratory or skin sensitization	No data available
Germ cell mutagenicity	No data available
Carcinogenicity	No data available
Reproductive toxicity	No data available
STOT-single exposure	No data available
STOT-repeated exposure	Kidneys and central nervous system
Aspiration hazard	No data available

Section 12 – Ecological Information

Environmental Toxicity	No Data Available
Aquatic Toxicity	No Data Available
Persistence and degradability	No Data Available
Bioaccumulative potential	No Data Available
Mobility in soil	Water soluble
Results of PBT and vPvT assessment	No data available

Section 13 – Disposal Considerations

Minimize waste as much as possible.

Disposal must be made according to state and federal regulations.

Section 14 – Transport Information

DOT (Ground)

Not regulated

IMDG

Not regulated

IATA

Not regulated

ADR

Not regulated

ADN

Not regulated

RID

Not regulated

Section 15 – Regulatory Information

15.1 Product related hazard information: The product has been classified and marked in accordance with regulations on hazardous materials.

Section 16 – Other information

Revision Date: June 3rd, 2014

To the best of our knowledge, the information contained herein is accurate. However, neither BioLegend, nor any of its subsidiaries assumes any liabilities whatsoever for the accuracy or completeness of the information contained herein. Final determination of suitability of any material is the sole responsibility of the user. All materials may present unknown hazards and should be used with caution. Although certain hazards are described herein, we cannot guarantee that these are the only hazards that exist.

Department issuing MSDS: Safety & Environment Department

Contact: Technical Service Representative