

IncuCyte™ Annexin V Red Reagent

1 IDENTIFICATION Includes product identifier; manufacturer or distributor name, address, phone number; emergency phone number; recommended use; restrictions on use.

Product code: Cat. No. 4641
 Product name: IncuCyte™ Annexin V Red Reagent
 Unit size: 1 vial, 100 tests

Manufacturer/supplier: Essen BioScience Inc.
 300 West Morgan Road
 Ann Arbor
 MI 48108
 USA

Emergency contact: +1 (734) 769-1600
 Fax contact: +1 (734) 769-7295

For research use only. Not intended for human or animal diagnostic or therapeutic uses.

2 HAZARD(S) IDENTIFICATION Includes all hazards regarding the chemical; required label elements.

GHS - Classification

Signal word: Not hazardous
 Health hazard: Not hazardous
 Physical hazards: Not hazardous

Hazard Statements

None

Precautionary Statements

None

Principle routes of exposure/Potential health effects

Eyes: Mild eye irritation in susceptible persons.
 Skin: Mild skin irritation.
 Inhalation: May be harmful by inhalation.
 Ingestion: Ingestion May be harmful if swallowed.

Specific effects

Carcinogenic effects: None.
 Mutagenic effects: None.
 Reproductive toxicity: None.
 Sensitization: No sensitization responses were observed.
 Target organ effects: No known effects under normal use conditions.
 HMIS: No information available.

3 COMPOSITION/INFORMATION ON INGREDIENTS Includes information on chemical ingredients; trade secret claims.

Chemical name	CAS-No	EINECS-No	Weight %
IncuCyte™ Annexin V Red Reagent	-	-	>99%

4 FIRST-AID MEASURES

Includes important symptoms/ effects, acute, delayed; required treatment.

- Skin contact: Wash off immediately with plenty of water. If symptoms occur, obtain medical advice.
 Eye contact: Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. Consult a physician if necessary.
 Ingestion: Never give anything by mouth to an unconscious person. If symptoms persist, call a physician. Do not induce vomiting without medical advice.
 Inhalation: Move to fresh air. If symptoms persist, call a physician. If not breathing, give artificial respiration.
 Notes to physician: Treat symptomatically.

5 FIRE-FIGHTING MEASURES

Lists suitable extinguishing techniques, equipment; chemical hazards from fire.

- Suitable extinguishing media: Water spray. Carbon dioxide (CO₂). Foam. Dry chemical.
 Special protective equipment for firefighters: Wear self-contained breathing apparatus and protective suit.

6 ACCIDENTAL RELEASE MEASURES

Lists emergency procedures; protective equipment; proper methods of containment and cleanup.

- Personal precautions: ELIMINATE all ignition sources (no smoking, flares, sparks or flames in immediate area). Use personal protection equipment. Avoid breathing vapors, mist or gas.
 Methods for cleaning up: Soak up with inert absorbent material. Sweep up and shovel into suitable containers for disposal. After cleaning, flush away traces with water.
 Environmental precautions: Prevent further leakage or spillage if safe to do so. Do not let product enter drains. See Section 12 for more information.

7 HANDLING AND STORAGE

Lists precautions for safe handling and storage, including incompatibilities.

- Handling: Avoid contact with skin and eyes. Always wear recommended Personal Protective Equipment. Avoid inhalation of vapor or mist. Avoid direct contact with substance.
 Storage: Keep away from heat, sparks, flame and other sources of ignition (i.e., pilot lights, electric motors and static electricity). Do not store near combustible materials. Keep in properly labeled containers.

**8 EXPOSURE CONTROLS/
PERSONAL PROTECTION**

Lists OSHA's Permissible Exposure Limits (PELs); ACGIH Threshold Limit Values (TLVs); and any other exposure limit used or recommended by the chemical manufacturer, importer, or employer preparing the SDS where available as well as appropriate engineering controls; personal protective equipment (PPE).

Chemical name	OSHA PEL	OSHA PEL (Ceiling)	ACGIH OEL (TWA)	ACGIH OEL (STEL)
IncuCyte™ Annexin V Red Reagent	None	None	None	None
Engineering measures:	Ensure adequate ventilation, especially in confined areas.			
Personal protective equipment:	Personal Protective Equipment requirements are dependent on the user institution's risk assessment and are specific to the risk assessment for each laboratory where this material may be used.			
Respiratory protection:	In case of insufficient ventilation wear suitable respiratory equipment.			
Hand protection:	Impervious gloves.			
Eye protection:	Safety glasses with side-shields.			
Skin and body protection:	Lightweight protective clothing.			
Hygiene measures:	Handle in accordance with good industrial hygiene and safety practice.			
Environmental exposure controls:	Prevent product from entering drains.			

9 PHYSICAL AND CHEMICAL PROPERTIES Lists the chemical's characteristics.

General information

Form: Solid.
 Appearance: No information available.
 Odor: No data available.
 Odor threshold: No data available.
 Boiling point/range: No data available.
 Melting point/range: No data available.
 Flash point: No data available.
 Autoignition temperature: No data available.
 Oxidizing properties: No information available.
 Water solubility: Miscible.
 Upper explosion limit: No data available.
 Lower explosion limit: No data available.
 Viscosity: No data available
 pH range: No data available.

10 STABILITY AND REACTIVITY Lists chemical stability and possibility of hazardous reactions.

Stability: Stable under normal conditions.
 Materials to avoid: No data available.
 Hazardous decomposition products: Hazardous decomposition products formed under fire conditions. Sulphur oxides. Carbon oxides.
 Polymerization: Hazardous polymerization does not occur.

11 TOXICOLOGICAL INFORMATION Includes routes of exposure; related symptoms, acute and chronic effects; numerical measures of toxicity

Acute toxicity

<u>Chemical name</u>	<u>LD50 (oral, rat/mouse)</u>	<u>LD50 (dermal, rat/rabbit)</u>	<u>LC50 (inhalation, rat/mouse)</u>
IncuCyte™ Annexin V Red Reagent	No data available	No data available	No data available

Principle routes of exposure/potential health effects

Eyes: Mild eye irritation in susceptible persons.
 Skin: Mild skin irritation in susceptible persons.
 Inhalation: No information available.
 Ingestion: No information available.
 Carcinogenic effects: None.
 Mutagenic effects: None.
 Reproductive toxicity: None.
 Sensitization: May cause sensitization in susceptible persons.

12 ECOLOGICAL INFORMATION List chemical's impact on the environment.

Ecotoxicity effects: No information available.
 Mobility: No information available.
 Biodegradation: No information available.
 Bioaccumulation: No information available.

13 DISPOSAL CONSIDERATIONS Prevent product from entering drains. Dispose of in accordance with local regulations.

14 TRANSPORT INFORMATION IATA.

Proper shipping name: No dangerous good in sense of these transport regulations.
 Hazard class: None.
 Subsidiary class: None.
 Packing group: None.
 UN-No: None.

15 REGULATORY INFORMATION Lists federal and state regulations.

Component	TSCA
IncuCyte™ Annexin V Red Reagent	-

US Federal Regulations

SARA 313: This product is not regulated by SARA.
 Clean air act, section 112 hazardous air pollutants (HAPs) (see 40 CFR 61): This product does not contain HAPs.

US State Regulations

California Proposition 65: This product does not contain chemicals listed under Proposition 65.
 WHMIS Hazard Class: Non-controlled.

Chemical name	Massachusetts RTK	New Jersey RTK	Pennsylvania RTK	Illinois RTK	Rhode Island RTK
IncuCyte™ Annexin V Red Reagent	-	-	-	-	-

California Proposition 65: This product does not contain any Proposition 65 chemicals.
 WHMIS Hazard Class: Non-controlled.



This product has been classified in accordance with the hazard criteria of the Controlled Products Regulations (CPR) and the MSDS contains all the information required by the CPR.

16 OTHER INFORMATION

Includes the date of preparation or last revision.

Date of preparation: March 3, 2016

Revision date:

Reason for revision:

For research use only. Not intended for human or animal diagnostic or therapeutic uses.

The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may present unknown hazards and should be used with caution. Since the Company cannot control the actual methods, volumes, or conditions of use, the Company shall not be held liable for any

damages or losses resulting from the handling or from contact with the product as described herein. THE INFORMATION IN THIS SDS DOES NOT CONSTITUTE A WARRANTY, EXPRESSED OR IMPLIED, INCLUDING ANY IMPLIED WARRANTY OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE.

END OF SAFETY DATA SHEET