

## IncuCyte® Cell Cycle Red/Green Lentivirus (EF-1 Alpha, Puro)

1. IDENTIFICATION	
Product code:	Catalog Number 4779
Product name:	IncuCyte® Cell Cycle Red/Green Lentivirus (EF-1 Alpha, Puro)
Unit size:	0.2 mL
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 IDENTIFICATION

Replication-defective lentiviral vectors, such as the 3rd generation vector provided in this product, are not known to cause any diseases in humans or animals. However, lentivirus particles still pose some biohazardous risk because they can transduce primary human cells and can integrate into the host cell genome thus posing some risk of insertional mutagenesis. For this reason, we highly recommend that you treat lentiviral stocks as Biosafety Level 2 (BSL-2,

BL-2) organisms and strictly follow all published BL-2 guidelines with proper waste decontamination. For more information about the BL-2 guidelines and lentivirus handling, we recommend referring to local documentation based on geography. The Essen BioScience 3rd generation HIV-based lentiviruses meet BL-2 requirements based on the criteria in the document, "Biosafety in Microbiological and Biomedical Laboratories", 5th Edition, published by the Centers for Disease Control (CDC).

This document may be downloaded at http://www.cdc.gov/biosafety/ publications/bmbl5/index.htm

## **Institutional Guidelines:**

Safety requirements for use and handling of lentiviruses may vary at individual institutions. We recommend consulting your institution's health and safety guidelines and/or officers prior to implementing the use of these reagents in your experiments.

Signal word:	Not hazardous.		
Hazard statement:	Not hazardous.		
Hazard pictogram:	None		
Principle routes of exposure/Potential health effects			
Eyes:	No information available.		
Skin:	No information available.		
Inhalation:	No information available.		
Ingestion:	May be harmful if swallowed.		
Specific effects			
Carcinogenic effects:	No information available.		
Mutagenic effects:	No information available.		
Reproductive toxicity:	No information available.		
Sensitization:	No information available.		
Target organ effects:	No information available.		
HMIS:	Health	Flammability	Reactivity
	0	0	0





3. COMPOSITION / INFORMATION ON INGREDIENTS		
	Describing at least in ince	
Hazardous /non-hazardous components	Recombinant lentivirus	
4. FIRST-AID MEASURES		
Skin contact:	Rinse with plenty of water. If symptoms arise, call a physician.	
Eye contact:	Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. If symptoms persist, call a physician.	
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		
Suitable extinguishing media:	Water spray, alcohol-resistant foam, dry chemical or CO <sub>2</sub> .	
Special protective equipment for fire fighters:	Wear self-contained breathing apparatus and protective suit.	
6. ACCIDENTAL RELEASE MEASURES		
Personal precautions:	Evacuate area. Wear self-contained breathing apparatus, rubber boots and heavy rubber gloves.	
Methods and materials for containment and cleaning up:	Absorb with sand or vermiculite, sweep up, place in a bag and hold for waste disposal. Avoid raising dust. Ventilate area and wash spill site after material pickup is complete.	
Environmental precautions:	Prevent product from entering drains.	
7. HANDLING AND STORAGE		
Handling:	Should be handled by trained personnel observing good laboratory practices.	
Storage:	Keep in properly labeled containers.	
8. EXPOSURE CONTROLS/PERSONAL PROTECTION		
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.	





9. PHYSICAL AND CHEMICAL PROPERTIES	
General information	
Form:	No information available.
Important health and safety environmental inf	ormation
Boiling point:	No information available.
Melting point/range:	No information available.
Flash point:	No information available.
Autoignition temperature:	No information available.
Oxidizing properties:	No information available.
Water solubility:	No data available.
10. STABILITY AND REACTIVITY	
Stability:	Stable under normal conditions.
Materials to avoid:	No dangerous reaction known under conditions of normal use.
Hazardous decomposition products:	No information available.
Polymerization:	Hazardous polymerization does not occur.
11. TOXICOLOGICAL INFORMATION	
Acute toxicity:	Not hazardous.
Principle routes of exposure/potential health ef	-
Eyes:	No information available.
Skin:	No information available.
Inhalation:	No information available.
Ingestion:	May be harmful by ingestion.
Carcinogenic effects:	No information available.
Mutagenic effects:	No information available.
Reproductive toxicity:	No information available.
Sensitization:	No information available.
Target organ effects:	No information available.
12. ECOLOGICAL INFORMATION	
Ecotoxicity effects:	No information available.
Mobility:	No information available.
Bioaccumulation:	Inherently biodegradable.
Bioaccumulation:	Does not bioaccumulate.
13. DISPOSAL CONSIDERATIONS	





14. TRANSPORTATION INFORMATION	
DOT (US)	
UN number:	3373
Hazard class:	6.2
Proper shipping name:	Biological substance, Category B
Reportable quantity (RQ):	No information available.
Poison inhalation hazard:	No
IMDG	
UN number:	3373
Hazard class:	6.2
Proper shipping name:	Biological substance, Category B
IATA	
UN number:	3373
Hazard class:	6.2
Proper shipping name:	Biological substance, Category B
15. REGULATORY INFORMATION	
US Federal Regulations	
SARA 313:	Not determined.
Clean air act, section 112 hazardous	Not listed.
air pollutants (HAPs) (see 40 CFR 61):	
US State Regulations	
California proposition 65:	This product does not contain chemicals listed under Proposition 65.
WHMIS hazard class:	Non-controlled.

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.

Date of preparation:		
ace of preparations	Date of preparation.	

## **END OF SAFETY DATA SHEET**

