

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

### SECTION 1: Identification of the substance/mixture and of the company/undertaking

#### 1.1 Product identifier

Trade name : Incucyte® Neurolight Orange Lentivirus

Product code : 4758

Substance name : Incucyte® Neurolight Orange Lentivirus

#### 1.2 Relevant identified uses of the substance or mixture and uses advised against

Recommended restrictions on use : Reserved for industrial and professional use., We recommend consulting your institution's health and safety guidelines and/or officers prior to implementing the use of these reagents in your experiments.

#### 1.3 Details of the supplier of the safety data sheet

Company : Sartorius Corporation  
565 Johnson Avenue  
11716 Bohemia

E-mail address of person responsible for the SDS : CA-FM Department: Environment@sartorius.com

#### 1.4 Emergency telephone number

UK/Europe/Global: +44.1865.407.333 (Sartorius 29003)  
United States: +1.866.928.0789 (Toll free), (Sartorius 29003)  
Canada: +1.800.579.7421 (Toll free), (Sartorius 29003)  
United States & Canada (geographic): +1.202.464.2554 (Sartorius 29003)  
Asia-Pacific/East & South East Asia (without China): +65.3158.1074 (Sartorius 29003)  
Japan: +81.3.4578.9341 (Sartorius 29003)  
Indonesia: 007.803.011.0293 (Toll free Indonesia), (Sartorius 29003)  
Malaysia: +60.3.6207.4347 (Sartorius 29003)  
Thailand: 001.800.120.666.751 (Toll free Thailand), (Sartorius 29003)  
India: 000.800.100.7479 (Toll free India) or +65.3158.1198 (Sartorius 29003)  
South Korea: +82.2.3479.8401 (Sartorius 29003)  
Australia: +61.2.8014.4558 (Sartorius 29003)  
New Zealand: +64.9.929.1483 (Sartorius 29003)  
Middle East/Africa: +44.1235.239671 (Sartorius 29003)

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

---

### SECTION 2: Hazards identification

#### 2.1 Classification of the substance or mixture

##### Classification (REGULATION (EC) No 1272/2008)

Not a hazardous substance or mixture according to the Globally Harmonised System (GHS).

#### 2.2 Label elements

##### Labelling REGULATION (EC) No 1272/2008

Not a hazardous substance or mixture according to the Globally Harmonised System (GHS).

Hazard pictograms :



Precautionary statements :

P261

Avoid breathing dust/ fume/ gas/ mist/ vapours/ spray.

P311

Call a POISON CENTER /doctor.

P272

Contaminated work clothing should not be allowed out of the workplace.

P281

Use personal protective equipment as required.

P280

Wear protective gloves/ eye protection/ face protection.

#### 2.3 Other hazards

We recommend consulting your institution's health and safety guidelines and/or officers prior to implementing the use of these reagents in your experiments.

Recombinant lentivirus/BSL-2 - BL-2

---

### SECTION 3: Composition/information on ingredients

#### 3.2 Mixtures

Chemical nature : Recombinant lentivirus/BSL-2 - BL-2

##### Hazardous components

Remarks : Recombinant lentivirus/BSL-2 - BL-2

---

## **Incucyte® Neurolight Orange Lentivirus**

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

---

### **SECTION 4: First aid measures**

#### **4.1 Description of first aid measures**

- |                         |   |
|-------------------------|---|
| General advice          | : No information available.   |
| If inhaled              | : If breathed in, move person into fresh air.<br>If symptoms persist, call a physician.   |
| In case of skin contact | : Wash off immediately with plenty of water.<br>If skin irritation persists, call a physician.  |
| In case of eye contact  | : Rinse immediately with plenty of lukewarm water, also under<br>the eyelids, for at least 15 minutes.<br>If eye irritation persists, consult a specialist. |
| If swallowed            | : Call a physician immediately.<br>DO NOT induce vomiting unless directed to do so by a physi-<br>cian or poison control center.                            |

#### **4.2 Most important symptoms and effects, both acute and delayed**

None known.

#### **4.3 Indication of any immediate medical attention and special treatment needed**

Treatment : No information available.

---

### **SECTION 5: Firefighting measures**

#### **5.1 Extinguishing media**

Suitable extinguishing media : Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.

#### **5.2 Special hazards arising from the substance or mixture**

Specific hazards during fire-fighting : No information available.

#### **5.3 Advice for firefighters**

Special protective equipment for firefighters : Wear self-contained breathing apparatus for firefighting if necessary.

Further information : The product itself does not burn.  
Use extinguishing measures that are appropriate to local circumstances and the surrounding environment.

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

---

### SECTION 6: Accidental release measures

#### 6.1 Personal precautions, protective equipment and emergency procedures

Personal precautions : Immediately evacuate personnel to safe areas.  
Use personal protective equipment.

#### 6.2 Environmental precautions

Environmental precautions : Do not let product enter drains.

#### 6.3 Methods and material for containment and cleaning up

Methods for cleaning up : Soak up with inert absorbent material (e.g. sand, silica gel, acid binder, universal binder, sawdust).

Biohazard decontamination/disinfection/inactivation.

#### 6.4 Reference to other sections

---

### SECTION 7: Handling and storage

#### 7.1 Precautions for safe handling

Advice on safe handling : Avoid formation of dust and aerosols.  
All processes must be supervised by specialists or authorised personnel.

Hygiene measures : Handle in accordance with good industrial hygiene and safety practice. Avoid contact with skin, eyes and clothing. Avoid contact with the skin and the eyes. Written instructions for handling must be available at the work place.

#### 7.2 Conditions for safe storage, including any incompatibilities

Requirements for storage areas and containers : Store in original container. Keep container tightly closed.

Storage class (TRGS 510) : 6.1D, Non-combustible, acute toxic Cat.3 / toxic hazardous materials or hazardous materials causing chronic effects

#### 7.3 Specific end use(s)

---

### SECTION 8: Exposure controls/personal protection

#### 8.1 Control parameters

#### 8.2 Exposure controls

Personal protective equipment

---

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

Eye protection	: Face-shield
Hand protection	
Remarks	: Rubber gloves
Skin and body protection	: Lab coat
	Rubber or plastic boots
Respiratory protection	: Self-contained closed-circuit breathing apparatus for escape
Protective measures	: Avoid contact with the skin and the eyes.
	When using do not eat or drink.

---

### SECTION 9: Physical and chemical properties

#### 9.1 Information on basic physical and chemical properties

Appearance	: liquid
Density	: Not applicable

#### 9.2 Other information

No data available

---

### SECTION 10: Stability and reactivity

#### 10.1 Reactivity

#### 10.2 Chemical stability

#### 10.3 Possibility of hazardous reactions

Hazardous reactions	: No dangerous reaction known under conditions of normal use. Stable under recommended storage conditions.
---------------------	---

#### 10.4 Conditions to avoid

Conditions to avoid	: Exposure to light.
---------------------	----------------------

#### 10.5 Incompatible materials

Materials to avoid	: Not applicable
--------------------	------------------

#### 10.6 Hazardous decomposition products

---

### SECTION 11: Toxicological information

#### 11.1 Information on toxicological effects

##### Further information

##### Product:

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

Remarks: biohazard  
Recombinant lentivirus/BSL-2 - BL-2

---

### SECTION 12: Ecological information

#### 12.1 Toxicity

No data available

#### 12.2 Persistence and degradability

No data available

#### 12.3 Bioaccumulative potential

No data available

#### 12.4 Mobility in soil

No data available

#### 12.5 Results of PBT and vPvB assessment

Not relevant

#### 12.6 Other adverse effects

**Product:**

Additional ecological information : Remarks: An environmental hazard cannot be excluded in the event of unprofessional handling or disposal.

---

### SECTION 13: Disposal considerations

#### 13.1 Waste treatment methods

Product : In accordance with local and national regulations.

---

### SECTION 14: Transport information

#### 14.1 UN number

ADR : UN 3373

IMDG : UN 3373

IATA : UN 3373

#### 14.2 UN proper shipping name

ADR : BIOLOGICAL SUBSTANCE, CATEGORY B

IMDG : BIOLOGICAL SUBSTANCE, CATEGORY B

IATA : BIOLOGICAL SUBSTANCE, CATEGORY B

---

---

## **Incucyte® Neurolight Orange Lentivirus**

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

### **14.3 Transport hazard class(es)**

<b>ADR</b>	: 6.2
<b>IMDG</b>	: 6.2
<b>IATA</b>	: 6.2

### **14.4 Packing group**

<b>ADR</b>	
Packing group	: Not assigned by regulation
Classification Code	: I4
Labels	: 6.2

<b>IMDG</b>	
Packing group	: II
Labels	: 6.2

<b>IATA</b>	
Packing group	: II
Labels	: 6.2

### **14.5 Environmental hazards**

<b>ADR</b>	
Environmentally hazardous	: no

<b>IMDG</b>	
Marine pollutant	: no

### **14.6 Special precautions for user**

Not applicable

### **14.7 Transport in bulk according to Annex II of MARPOL 73/78 and the IBC Code**

Not applicable for product as supplied.

---

## **SECTION 15: Regulatory information**

### **15.1 Safety, health and environmental regulations/legislation specific for the substance or mixture**

Water contaminating class (Germany)	: WGK 3 highly hazardous to water
-------------------------------------	-----------------------------------

Other regulations	: Restricted to professional users. Attention - Avoid exposure - obtain special instructions before use.
-------------------	--

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## Incucyte® Neurolight Orange Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: -
1.0	16.09.2021	SAG-R0002651	Date of first issue: 16.09.2021

---

### 15.2 Chemical safety assessment

---

## SECTION 16: Other information

### Further information

Training advice : Provide adequate information, instruction and training for operators.

The information provided in this Safety Data Sheet is correct to the best of our knowledge, information and belief at the date of its publication. The information given is designed only as a guidance for safe handling, use, processing, storage, transportation, disposal and release and is not to be considered a warranty or quality specification. The information relates only to the specific material designated and may not be valid for such material used in combination with any other materials or in any process, unless specified in the text.

DE / EN