

SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

Incucyte® Neurolight Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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

1.1 Product identifier

Trade name : Incucyte® Neurolight Red Lentivirus

Product code : 4584

Substance name : Incucyte® Neurolight Red 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 Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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 Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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.
If symptoms persist, call a physician. |
| In case of skin contact | : | Wash off immediately with plenty of water.
If skin irritation persists, call a physician. |
| In case of eye contact | : | Rinse immediately with plenty of lukewarm water, also under
the eyelids, for at least 15 minutes.
If eye irritation persists, consult a specialist. |
| If swallowed | : | Call a physician immediately.
DO NOT induce vomiting unless directed to do so by a physi-
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 car- bon 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 nec- essary.
Further information	:	The product itself does not burn. Use extinguishing measures that are appropriate to local cir- cumstances and the surrounding environment.

SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

Incucyte® Neurolight Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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

Incucyte® Neurolight Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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 Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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 Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

14.3 Transport hazard class(es)

ADR	: 6.2
IMDG	: 6.2
IATA	: 6.2

14.4 Packing group

ADR	
Packing group	: Not assigned by regulation
Classification Code	: I4
Labels	: 6.2

IMDG	
Packing group	: II
Labels	: 6.2

IATA	
Packing group	: II
Labels	: 6.2

14.5 Environmental hazards

ADR	
Environmentally hazardous	: no

IMDG	
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 hazard 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 Red Lentivirus

Version	Revision Date:	SDS Number:	Date of last issue: 18.08.2021
1.2	09.02.2022	SAG-R0001869	Date of first issue: 30.03.2020

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