

SAFETY DATA SHEET

CellBrite Cytoplasmic Membrane Dyes

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

1.1. Product identifier

Trade name

CellBrite Cytoplasmic Membrane Dyes

Other names / Synonyms

30021: CellBrite® Green Cytoplasmic Membrane Dye, 1 mL 30022: CellBrite® Orange Cytoplasmic Membrane Dye, 1 mL 30023: CellBrite® Red Cytoplasmic Membrane Dye, 1 mL Product no.

30021, 30022, 30023

Other means of identification Index No.: 603-002-00-5 EC No.: 200-578-6

CAS No.: 64-17-5

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

Relevant identified uses of the substance or mixture

Use as laboratory reagent. For research use only. Not for food, drug, household, or cosmetic use. Restricted to professional users.

Use descriptors (REACH)

Sectors of use	Description
SU 24	Scientific research and development
Product category	Description
PC 21	Laboratory Chemicals
Process category	Description
PROC 15	Use as laboratory reagent
ses advised against	

None known.

1.3. Details of the supplier of the safety data sheet

Company and address

```
Biotium, Inc.
```

46117 Landing Parkway CA 94538 Fremont USA T: +1 510-265-1027 Fax: +1 510-265-1352 http://www.biotium.com E-mail techsupport@biotium.com Revision 02/05/2024 SDS Version 1.0 Date of previous version 02/05/2024 (1.0) 1.4. Emergency telephone number

Anti-poison center: +32 (0) 70 245 245 (every day, 24 hours a day) See also section 4 "First aid measures".



SECTION 2: Hazards identification

```
Classified according to Regulation (EC) No. 1272/2008 (CLP).
2.1. Classification of the substance or mixture
  Flam. Liq. 2; H225, Highly flammable liquid and vapour.
2.2. Label elements
  Hazard pictogram(s)
  Signal word
      Danger
  Hazard statement(s)
     Highly flammable liquid and vapour. (H225)
  Precautionary statement(s)
     General
      Prevention
        Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking. (P210)
        Keep container tightly closed. (P233)
      Response
       In case of fire: Use water mist/carbon dioxide/alcohol-resistant foam to extinguish. (P370+P378)
      Storage
        Store in a well-ventilated place. Keep cool. (P403+P235)
      Disposal
        Dispose of contents/container in accordance with local regulation
        (P501)
  Hazardous substances
     ethanol; ethyl alcohol
  Additional labelling
     Not applicable.
2.3. Other hazards
  Additional warnings
```

This mixture/product does not contain any substances known to fulfil the criteria for PBT and vPvB classification. This product does not contain any substances considered to be endocrine disruptors in accordance with the criteria set out in Commission Delegated Regulation (EU) 2017/2100 or Commission Regulation (EU) 2018/605.

SECTION 3: Composition/information on ingredients

3.1. ▼Substances

Product/substance	Identifiers	% w/w	Classification	Note
ethanol;ethyl alcohol	CAS No.: 64-17-5 EC No.: 200-578-6 REACH: Index No.: 603-002-00-5	>99%%	Flam. Liq. 2, H225	

3.2. Mixtures

Not applicable. This product is a substance.

See full text of H-phrases in section 16. Occupational exposure limits are listed in section 8, if these are available.

Other information



SECTION 4: First aid measures

4.1. Description of first aid measures

General information

In the case of accident: Contact a doctor or casualty department – bring the label or this safety data sheet. The doctor can contact the Anti-poison center: +32 (0) 70 245 245 (every day, 24 hours a day)

Contact a doctor if in doubt about the injured person's condition or if the symptoms persist. Never give an unconscious person water or other drink.

Inhalation

Upon breathing difficulties or irritation of the respiratory tract: Bring the person into fresh air and stay with him/her.

Skin contact

Remove contaminated clothing and shoes immediately. Ensure to wash exposed skin thoroughly with water and soap. Skin cleanser can be used. DO NOT use solvents or thinners.

Eye contact

If in eyes: Flush eyes with water or saline water (20-30 °C) for at least 5 minutes. Remove contact lenses. Seek medical assistance and continue flushing during transport.

Ingestion

If the person is conscious, rinse the mouth with water and stay with the person. Never give the person anything to drink.

In case of malaise, seek medical advice immediately and bring the safety data sheet or label from the product. Do not induce vomiting, unless recommended by the doctor. Have the person lean forward with head down to avoid inhalation of or choking on vomited material.

Burns

Rinse with water until pain stops then continue to rinse for 30 minutes.

4.2. Most important symptoms and effects, both acute and delayed

Neurotoxic effects: This product contains organic solvents, which may cause adverse effects to the nervous system. Symptoms of neurotoxicity include: loss of appetite, headache, dizziness, ringing in ears, tingling sensations of skin, sensitivity to the cold, cramps, difficulty in concentrating, tiredness, etc. Repeated exposure to solvents can result in the breaking down of the skin's natural fat layer and may result in an increased absorption potential of other hazardous substances at the area of exposure.

4.3. Indication of any immediate medical attention and special treatment needed

Treat symptomatically.

Information to medics

Bring this safety data sheet or the label from this product.

SECTION 5: Firefighting measures

5.1. Extinguishing media

Suitable extinguishing media: Alcohol-resistant foam, carbon dioxide, powder, water mist. Unsuitable extinguishing media: Waterjets should not be used, since they can spread the fire.

5.2. Special hazards arising from the substance or mixture

Highly flammable liquid and vapour.

In use may form flammable/explosive vapour-air mixture.

Fire will result in dense smoke. Exposure to combustion products may harm your health. Closed containers, which are exposed to fire, should be cooled with water. Do not allow fire-extinguishing water to enter the sewage system and nearby surface waters.

If the product is exposed to high temperatures, e.g. in the event of fire, dangerous decomposition compounds are produced. These are:

Carbon oxides (CO / CO2)

5.3. Advice for firefighters

Wear self-contained breathing apparatus and protective clothing to prevent contact. Upon direct exposure contact the anti-poison center: +32 (0) 70 245 245 (every day, 24 hours a day) in order to obtain further advice.

SECTION 6: Accidental release measures

6.1. Personal precautions, protective equipment and emergency procedures

Storages not yet ignited must be cooled by water mist. Remove flammable materials if conditions allow it. Ensure



sufficient ventilation. Ensure adequate ventilation, especially in confined areas. Contaminated areas may be slippery.

6.2. Environmental precautions

Avoid discharge to lakes, streams, sewers, etc. Keep unauthorized persons away from the spill

Reep unautionized persons away from the spin

6.3. Methods and material for containment and cleaning up

Contain and collect spillage with non-combustible, absorbent material e.g. sand, earth, vermiculite or diatomaceous earth and place in container for disposal according to local regulations. Wherever possible cleaning should be performed with normal cleaning agents. Avoid use of solvents.

6.4. Reference to other sections

See section 13 "Disposal considerations" on handling of waste.

See section 8 "Exposure controls/personal protection" for protective measures.

SECTION 7: Handling and storage

7.1. Precautions for safe handling

Ground and bond container and receiving equipment. Use explosion-proof [electrical/lighting/ventilating] equipment. Use non-sparking tools. Take action to prevent static discharges. Avoid contact during pregnancy and while nursing. Smoking, drinking and consumption of food is not allowed in the work area. See section 8 "Exposure controls/personal protection" for information on personal protection.

7.2. Conditions for safe storage, including any incompatibilities

Containers that have been opened must be carefully resealed and kept upright to prevent leakage.

Take action to prevent static discharges.

Must be stored in a cool and well-ventilated area, away from possible sources of ignition.

Recommended storage material

Always store in containers of the same material as the original container.

Storage temperature

Keep container tightly closed in a dry and well-ventilated place. Refrigerator 2°C to 8°C. Protect from light.

Incompatible materials

No specific requirements

7.3. Specific end use(s)

This product should only be used for applications quoted in section 1.2.

SECTION 8: Exposure controls/personal protection

8.1. Control parameters

ethanol;ethyl alcohol Long term exposure limit (8 hours) (mg/m³): 1907 Long term exposure limit (8 hours) (ppm): 1000

List of the limit values for exposure to chemical agents (KB van May 2021).

DNEL

Duration: Long term – Systemic effects - Workers	Route of exposure:	DNEL: 343 mg/kg bw/day
		5 5 ,
Long term – Systemic effects - Workers	Inhalation	380 mg/m ³
Short term – Local effects - Workers	Inhalation	1900 mg/m³
PNFC		
ethanol;ethyl alcohol		
Route of exposure:	Duration of Exposure:	PNEC:



Freshwater	960 µg/L
Freshwater sediment	3.6 mg/kg
Intermittent release (freshwater)	2.75 mg/L
Marine water	790 μg/L
Marine water sediment	2.9 mg/kg
Predators	380-720 mg/kg
Sewage treatment plant	580 mg/L
Soil	630 µg/kg

8.2. Exposure controls

Compliance with the given occupational exposure limits values should be controlled on a regular basis.

General recommendations

Smoking, drinking and consumption of food is not allowed in the work area.

Exposure scenarios

There are no exposure scenarios implemented for this product.

Exposure limits

Professional users are subjected to the legally set maximum concentrations for occupational exposure. See occupational hygiene limit values above.

Appropriate technical measures

The formation of vapours must be kept at a minimum and below current limit values (see above). Installation of a local exhaust system if normal air flow in the work room is not sufficient is recommended. Ensure eyewash and emergency showers are clearly marked.

Apply standard precautions during use of the product. Avoid inhalation of vapours.

Hygiene measures

In between use of the product and at the end of the working day all exposed areas of the body must be washed thoroughly. Pay special attention to hands, forearms and face.

Measures to avoid environmental exposure

No specific requirements.

Individual protection measures, such as personal protective equipment

Generally

Use only CE marked protective equipment.

Respiratory Equipment

Туре	Class	Colour	Standards	
Not required; ex case of aerosol formation.	xcept in		N/A	

Skin protection

Recommended	Type/Category	Standards	
No specific requirements.	-	-	
Hand protection			

Material	Glove thickness (mm)	Breakthrough time (min.)	Standards	
Handle with gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory	2			



Material	Glove thickness (mm)	Breakthrough time (min.)	Standards
practices. Wash and dry hands.			
ye protection			
Туре	Standards		
Safety glasses with side-shields conforming to EN166. Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166(EU).	EN166		
CTION 9: Physical and ch	emical properties		
. Information on basic pł Physical state	nysical and chemical pro	operties	
Liquid			
Colour Testing not relevant o	r not possible due to th	e nature of the product.	
Odour / Odour threshold Testing not relevant o		e nature of the product.	
рН			
l esting not relevant o Density (g/cm³)	r not possible due to th	e nature of the product.	
Testing not relevant o	r not possible due to th	e nature of the product.	
Kinematic viscosity Testing not relevant o	r not possible due to th	e nature of the product.	
Particle characteristics Does not apply to liqu	ide		
ase changes	ius.		
Melting point/Freezing p		a mature of the preduct	
Softening point/range (w		e nature of the product.	
Does not apply to liqu			
Boiling point (°C) Testing not relevant o	r not possible due to th	e nature of the product.	
Vapour pressure			
Relative vapour density	ir not possible due to th	e nature of the product.	
Testing not relevant o		e nature of the product.	
Decomposition temperat Testing not relevant o		e nature of the product.	
ta on fire and explosion h			
Flash point (°C) Testing not relevant o	r not possible due to th	e nature of the product.	
Flammability (°C)			
The material is ignitat Auto-ignition temperatur			
Testing not relevant o	r not possible due to th	e nature of the product.	
Lower and upper explosi Testing not relevant o		e nature of the product.	
ubility		P	
Solubility in water			



n-octanol/water coefficient (LogKow) Testing not relevant or not possible due to the nature of the product. Solubility in fat (g/L) Testing not relevant or not possible due to the nature of the product. 9.2. Other information Other physical and chemical parameters No data available. Oxidizing properties Testing not relevant or not possible due to the nature of the product. SECTION 10: Stability and reactivity 10.1. Reactivity No data available. 10.2. Chemical stability The product is stable under the conditions, noted in section 7 "Handling and storage". 10.3. Possibility of hazardous reactions None known. 10.4. Conditions to avoid Do not expose to any forms of heat (e.g. solar radiation). May lead to excess pressure. Heat, flames and sparks. 10.5. Incompatible materials No specific requirements 10.6. Hazardous decomposition products The product is not degraded when used as specified in section 1. SECTION 11: Toxicological information 11.1. Information on hazard classes as defined in Regulation (EC) No 1272/2008 Acute toxicity Based on available data, the classification criteria are not met. Skin corrosion/irritation Based on available data, the classification criteria are not met. Serious eye damage/irritation Based on available data, the classification criteria are not met. **Respiratory sensitisation** Based on available data, the classification criteria are not met. Skin sensitisation Based on available data, the classification criteria are not met. Germ cell mutagenicity Based on available data, the classification criteria are not met. Carcinogenicity Based on available data, the classification criteria are not met. Reproductive toxicity Based on available data, the classification criteria are not met. STOT-single exposure Based on available data, the classification criteria are not met. STOT-repeated exposure Based on available data, the classification criteria are not met. Aspiration hazard Based on available data, the classification criteria are not met. 11.2. Information on other hazards Long term effects Neurotoxic effects: This product contains organic solvents, which may cause adverse effects to the nervous system. Symptoms of neurotoxicity include: loss of appetite, headache, dizziness, ringing in ears, tingling sensations of skin, sensitivity to the cold, cramps, difficulty in concentrating, tiredness, etc. Repeated exposure to solvents can result in the breaking down of the skin's natural fat layer and may result in an increased absorption potential of other hazardous substances at the area of exposure.



Endocrine disrupting properties

This mixture/product does not contain any substances known to have hormone-disrupting properties in relation to health.

Other information

None known.

SECTION 12: Ecological information

12.1. Toxicity

No data available.

12.2. Persistence and degradability

Based on available data, the classification criteria are not met.

12.3. Bioaccumulative potential

Based on available data, the classification criteria are not met.

- 12.4. Mobility in soil
 - No data available.
- 12.5. Results of PBT and vPvB assessment

This mixture/product does not contain any substances known to fulfil the criteria for PBT and vPvB classification.

12.6. Endocrine disrupting properties

This mixture/product does not contain any substances considered to have endocrine-disrupting properties in relation to the environment.

12.7. Other adverse effects None known.

SECTION 13: Disposal considerations

13.1. Waste treatment methods

Product is covered by the regulations on hazardous waste. HP 3 - Flammable Dispose of contents/container to an approved waste disposal plant. Commission Regulation (EU) No 1357/2014 of 18 December 2014 on waste. EWC code Not applicable.

Contaminated packing

Packaging containing residues of the product must be disposed of similarly to the product.

SECTION 14: Transport information

	14.1 UN / ID	14.2 UN proper shipping name	14.3 Hazard class(es)	14.4 PG*	14.5 Env**	Other information:
ADR	UN1170	ETHANOL SOLUTION (ETHYL ALCOHOL SOLUTION)	Transport hazard class: 3 Label: 3 Classification code: F1	Π	No	Limited quantities: 1 L Tunnel restriction code: (D/E) See below for additional information.
IMDG	UN1170	ETHANOL SOLUTION (ETHYL ALCOHOL SOLUTION)	Transport hazard class: 3 Label: 3 Classification code: F1	Π	No	Limited quantities: 1 L EmS: F-E S-D See below for additional information.
IATA	UN1170	ETHANOL SOLUTION (ETHYL ALCOHOL	Transport hazard class: 3	II	No	See below for



14.1 14.2	14.3	14.4	14.5	Other
UN / ID UN proper shipping name	Hazard class(es)	PG*	Env**	information:
SOLUTION)	Label: 3 Classification code: F1			additional information.

* Packing group

** Environmental hazards

Additional information

ADR / See Table A, section 3.2.1 for any information on special provisions, requirements, or warnings in connection with transport. See section 5.4.3, for instructions in writing regarding mitigation of damages in relation to incidents or accidents during transport.

IMDG / See section 3.2.1, for any information on special provisions, requirements, or warnings in connection with transport.

IATA / See Table 4.2 for any information on special provisions, requirements, or warnings in connection with transport.

This product is within scope of the regulations of transport of dangerous goods.

14.6. Special precautions for user

Not applicable.

14.7. Maritime transport in bulk according to IMO instruments No data available.

SECTION 15: Regulatory information

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

Restrictions for application

Restricted to professional users.

People under the age of 18 shall not be exposed to this product.

Pregnant women and women breastfeeding must not be exposed to this product. The risk, and possible technical precautions or design of the workplace needed to eliminate exposure, must be considered.

Demands for specific education

No specific requirements.

SEVESO - Categories / dangerous substances

P5c - FLAMMABLE LIQUIDS, Qualifying quantity (lower-tier): 5.000 tonnes / (upper-tier): 50.000 tonnes

REACH, Annex XVII

ethanol; ethyl alcohol is subject to REACH restrictions, REACH annex XVII (entry 40).

Additional information

Not applicable.

Sources

Labor Law of 24-01-1985, updated 31-12-2020.

5 June 2015 - Cooperation agreement between the Federal State, the Flemish Region, the Walloon Region and the Brussels-Capital Region concerning the control of hazards related to major accidents involving dangerous substances.

Commission Regulation (EU) No 1357/2014 of 18 December 2014 on waste.

Regulation (EC) No 1272/2008 of the European Parliament and of the Council of 16 December 2008 on

classification, labelling and packaging of substances and mixtures (CLP).

Regulation (EC) No 1907/2006 of the European Parliament and of the Council of 18 December 2006 concerning the Registration, Evaluation, Authorisation and Restriction of Chemicals (REACH).

15.2. Chemical safety assessment

No

SECTION 16: Other information

Full text of H-phrases as mentioned in section 3

H225, Highly flammable liquid and vapour.

The full text of identified uses as mentioned in section 1



SU 24 = Scientific research and development PROC 15 = Use as laboratory reagent PC 21 = Laboratory Chemicals Abbreviations and acronyms ADN = European Provisions concerning the International Carriage of Dangerous Goods by Inland Waterway ADR = The European Agreement concerning the International Carriage of Dangerous Goods by Road ATE = Acute Toxicity Estimate BCF = Bioconcentration Factor CAS = Chemical Abstracts Service CE = Conformité Européenne (European conformity) CLP = Classification, Labelling and Packaging Regulation [Regulation (EC) No. 1272/2008] CSA = Chemical Safety Assessment CSR = Chemical Safety Report DMEL = Derived Minimal Effect Level DNEL = Derived No Effect Level EINECS = European Inventory of Existing Commercial chemical Substances ES = Exposure Scenario EUH statement = CLP-specific Hazard statement EuPCS = European Product Categorisation System EWC = European Waste Catalogue GHS = Globally Harmonized System of Classification and Labelling of Chemicals IARC = International Agency for Research on Cancer (IARC) IATA = International Air Transport Association IBC = Intermediate Bulk Container IMDG = International Maritime Dangerous Goods LogPow = logarithm of the octanol/water partition coefficient MARPOL = International Convention for the Prevention of Pollution From Ships, 1973 as modified by the Protocol of 1978. ("Marpol" = marine pollution) OECD = Organisation for Economic Co-operation and Development PBT = Persistent, Bioaccumulative and Toxic PNEC = Predicted No Effect Concentration RID = The Regulations concerning the International Carriage of Dangerous Goods by Rail RRN = REACH Registration Number SCL = A specific concentration limit SVHC = Substances of Very High Concern STOT-RE = Specific Target Organ Toxicity - Repeated Exposure STOT-SE = Specific Target Organ Toxicity - Single Exposure TWA = Time weighted average UN = United Nations UVBC = Unknown or variable composition, complex reaction products or of biological materials VOC = Volatile Organic Compound vPvB = Very Persistent and Very Bioaccumulative Additional information The classification of the mixture in regard to physical hazards has been based on experimental data. The information provided above is believed to be correct to our best knowledge, but does not purport to be all

inclusive, and shall be used only as a quide. This material is sold for research purposes only and is not required to appear on the TSCA inventory. It is not intended for food, drug, household, agricultural or cosmetic use. Its use must be supervised by a technically qualified individual experienced in handling potentially hazardous chemicals. Biotium shall not be held liable for any damage resulting from handling or contact with the above product.

The safety data sheet is validated by

Eric Torres

Other

A change (in proportion to the last essential change (first cipher in SDS version, see section 1)) is marked with a triangle.

The information in this safety data sheet applies only to this specific product (mentioned in section 1) and is not necessarily correct for use with other chemicals/products.

It is recommended to hand over this safety data sheet to the actual user of the product. Information in this safety data sheet cannot be used as a product specification.

Country-language: BE-en