

SAFETY DATA SHEET

CellBrite Cytoplasmic Membrane Dyes

SECTION 1: Identification

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 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. Uses 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 SDS date 5/2/2024 **SDS Version** 1.0 Date of previous version 5/2/2024 (1.0) 1.4. Emergency telephone number Contact the poison control at 1-800-222-1222 (24/7) or use the webPOISONCONTROL® (triage.webpoisoncontrol.org) to get specific guidance for your case See also section 4 "First aid measures". SECTION 2: Hazard(s) identification **OSHA/HCS** status This material is considered hazardous by the OSHA Hazard Communication Standard (29 CFR 1910.1200) 2.1. Classification of the substance or mixture Flam. Lig. 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) |
| 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. |
| SECTION 3: Composition/Information on Ingredients |
| |

3.1. ▼Substances

| Product/substance | Identifiers | % w/w | Classification | Note |
|-----------------------|------------------|-------|--------------------|------|
| ethanol;ethyl alcohol | CAS No.: 64-17-5 | >99%% | Flam. Liq. 2, H225 | |

3.2. Mixtures

Not applicable. This product is a substance.

Where the concentration of an ingredient is expressed as a range the exact concentration has been withheld as a trade secret.

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

If breathing is irregular, drowsiness, loss of consciousness or cramps: Call 911 and give immediate treatment (first aid).

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: Fire-fighting 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 Poison Help Line on 1-800-222-1222 (24/7) 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

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 Short term exposure limit (STEL) (ACGIH TLV) (ppm): 1000 Long term exposure limit (OSHA Table Z-1) (mg/m³): 1900 Long term exposure limit (OSHA Table Z-1) (ppm): 1000

Part 1910 - Occupational Safety and Health Standards (29 CFR 1910.1000 TABLE Z-1 - Limits for Air Contaminants)

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 protective equipment with a recognized certification mark, e.g. the UL mark. Respiratory Equipment



| Туре | Class | Colour | Standards | |
|---|---|-----------------------------|-----------|--|
| Not required; except in case of aerosol formation. | | | N/A | |
| Skin protection | | | | |
| Recommended | Type/Category | Standar | ds | |
| 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 practices. Wash and dry hands. | | | | |
| Eye 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 | | | |
| ECTION 9: Physical and ch | nemical properties | | | |
| Odour Testing not relevant o Odour threshold (ppm) | nysical and chemical pro or not possible due to th or not possible due to th or not possible due to th | ne nature of the produc | :t. | |
| pH Testing not relevant o | or not possible due to th | ne nature of the produc | ct. | |
| Density (g/cm³) | or not possible due to th | | | |



Phase changes Melting point (°F) Testing not relevant or not possible due to the nature of the product. Softening point/range (waxes and pastes) (°F) Does not apply to liquids. Boiling point (°F) Testing not relevant or not possible due to the nature of the product. Vapour pressure Testing not relevant or not possible due to the nature of the product. Relative vapour density Testing not relevant or not possible due to the nature of the product. Decomposition temperature (°F) Testing not relevant or not possible due to the nature of the product. Data on fire and explosion hazards Flash point (°F) Testing not relevant or not possible due to the nature of the product. Flammability (°F) The material is ignitable. Auto-ignition temperature (°F) Testing not relevant or not possible due to the nature of the product. Explosion limits (% v/v) Testing not relevant or not possible due to the nature of the product. Solubility Solubility in water Soluble n-octanol/water coefficient (LogKow) Testing not relevant or not possible due to the nature of the product. Solubility in fat (q/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 toxicological effects

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.

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.

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. Other adverse effects None known.

SECTION 13: Disposal considerations

RCRA Hazardous waste ("P" and "U" list) (40 CFR 261)

None of the components are listed

Specific labelling

Contaminated packing

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

SECTION 14: Transport information

| | 14.1 14.2 | 14.3 | 14.4 | 14.5 | Other |
|-----|---|--|------|-------|--------------------------------------|
| | UN / ID UN proper shipping name | Hazard class(es) | PG* | Env** | information: |
| DOT | UN1170 ETHANOL SOLUTION (ETHYL ALCOHOL SOLUTION) | Transport hazard class: 3 Label: 3 Classification code: F1 | Π | No | Limited quantities: 1 L Tunnel |



| | 14.1 | 14.2 | 14.3 | 14.4 | 14.5 | Other |
|------|---------|--|--|------|-------|---|
| | UN / ID | UN proper shipping name | Hazard class(es) | PG* | Env** | information: |
| | | | | | | 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. |
| ΙΑΤΑ | UN1170 | ETHANOL SOLUTION (ETHYL ALCOHOL SOLUTION) | Transport hazard class: 3 Label: 3 Classification code: F1 | Π | No | See below for additional information. |

* Packing group

** Environmental hazards

Additional information

DOT / See § 172.101 Hazardous Materials Table for any information on special provisions, requirements, or warnings in connection with transport. See § 172.602, 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. Transport in bulk according to Annex II of Marpol and the IBC Code No data available.

SECTION 15: Regulatory information

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

15.2. U.S. Federal regulations

TSCA (the non-confidential portion) ethanol;ethyl alcohol is listed

Clean Air Act

None of the components are listed

EPCRA Section 302

None of the components are listed

EPCRA Section 304

None of the components are listed

EPCRA section 313

None of the components are listed

CERCLA

None of the components are listed

State regulations

California / Prop. 65

None of the components are listed Massachusetts / Right To Know Act



ethanol; ethyl alcohol is listed New Jersey / Right To Know Act ethanol; ethyl alcohol / Substance number: 0844 ethanol; ethyl alcohol is on the Special Health Hazard Substance List New York / Right To Know Act ethanol; ethyl alcohol is listed ethanol; ethyl alcohol is regulated with a Treshold Reporting Quantity (TRQ) of: 0 pounds Pennsylvania / Right To Know Act ethanol; ethyl alcohol is listed 15.4. Restrictions for application Restricted to professional users. 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. 15.5. Demands for specific education No specific requirements. 15.6. Additional information Not applicable. 15.7. Chemical safety assessment No 15.8. Sources OSHA Hazard Communication Standard (29 CFR 1910.1200) **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 None known. Abbreviations and acronyms ACGIH = American Conference of Governmental Industrial Hygienists 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 CERCLA = Comprehensive Environmental Response Compensation and Liability Act DOT = Department of Transportation EINECS = European Inventory of Existing Commercial chemical Substances EPCRA = Emergency Planning and Community Right-To-Know Act GHS = Globally Harmonized System of Classification and Labelling of Chemicals HCIS = Hazardous Chemical Information System HNOC = Hazards Not Otherwise Classified IARC = International Agency for Research on Cancer IATA = International Air Transport Association 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) NFPA = National Fire Protection Association NIOSH = National Institute for Occupational Safety and Health OECD = Organisation for Economic Co-operation and Development OSHA = Occupational Safety and Health Administration PBT = Persistent, Bioaccumulative and Toxic RCRA = Resource Conservation and Recovery Act RID = The Regulations concerning the International Carriage of Dangerous Goods by Rail RRN = REACH Registration Number SARA = Superfund Amendments and Reauthorization Act SCL = A specific concentration limit.



STEL = Short-term exposure limits

STOT-RE = Specific Target Organ Toxicity - Repeated Exposure

STOT-SE = Specific Target Organ Toxicity - Single Exposure

TSCA = The Toxic Substances Control Act

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 guide. 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: US-en